

SDMA ELISA

Manufactured by Immundiagnostik AG.

ALX-850-331-KI01

(Version: 2019-01-02)

Global Headquarters

ENZO LIFE SCIENCES, INC.

10 Executive Blvd.
Farmingdale, NY 11735
Ph: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

European Sales Office

ENZO LIFE SCIENCES (ELS) AG

Industriestrasse 17
CH-4415 Lausen, Switzerland
Ph: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com

LOCAL EUROPEAN OFFICES

Belgium, The Netherlands & Luxembourg

Enzo Life Sciences BVBA
Avenue Louise 65 / Box 11
1050 Bruxelles
Belgium
Ph: +32 3 466 0420
Fax: +32 3 808 7033
info-be@enzolifesciences.com

Germany

Enzo Life Sciences GmbH
Basler Strasse 57a
DE-79540 Lörrach
Germany
Ph: +49 7621 5500 526
Fax: +49 7621 5500 527
info-de@enzolifesciences.com

France

Enzo Life Sciences (ELS) AG
Branch Office Lyon
13, avenue Albert Einstein,
F-69100 Villeurbanne, France
Ph: +33 472 440 655
Fax: +33 481 680 254
info-fr@enzolifesciences.com

UK & Ireland

Enzo Life Sciences (UK) Ltd.
1 Colleton Crescent
Exeter EX2 4DG
Ph: 0845 601 1488 (UK customers)
Ph: +44 1392 825900
Fax: +44 1392 825910
info-uk@enzolifesciences.com

www.enzolifesciences.com





For laboratory use only. Not for human or diagnostic use.

Manual

SDMA ELISA

For the in vitro determination of SDMA in human, canine and feline serum, EDTA and Li-heparin plasma

Valid from 2019-01-02

REF **K 7780**  **96**  **+2°C**  **+8°C**  **-20°C**

STD
CTRL 1
CTRL 2

IVD



Immundiagnostik AG, Stubenwald-Allee 8a, 64625 Bensheim, Germany

Tel.: +49 6251 70190-0

Fax: + 49 6251 849430

e.mail: info@immundiagnostik.com

www.Immundiagnostik.com

1. INTENDED USE

This Immundiagnostik assay is intended for the quantitative determination of SDMA in human, canine and feline serum, EDTA and Li-heparin plasma. For *in vitro* diagnostic use only.

2. INTRODUCTION

The dosage of most drugs must be adapted in renal insufficiency, making accurate assessment of renal function a prerequisite in clinical medicine. Furthermore, even a modest decline in renal function has been recognized as a cardiovascular risk.

In clinical practice serum creatinine is typically used to assess renal function, but this serum creatinine does not increase at modest decline in renal function. Consequently, there is an ongoing search for suitable endogenous markers of renal function.

SDMA is a methylated derivative of L-arginine which is strictly eliminated by renal extraction, thus the SDMA plasma level is strongly correlated to renal function. In 18 studies with more than 2136 patients systemic SDMA concentrations correlated highly with inulin clearance, as well as with various clearance estimates combined and serum creatinine. With respect to this, SDMA exhibits properties of a reliable marker of renal dysfunction.

Moreover, there are hints that increased SDMA correlates with total sequential organ failure, indicating both renal and hepatic failure, and with an increased cardiovascular risk.

Indication

- Renal failure
- Cardiovascular risk in renal dysfunction
- Hypertension in renal dysfunction

3. MATERIAL SUPPLIED

Cat. No.	Label	Kit Components	Quantity
K 7780	PLATE	Microtiter plate, pre-coated	12 x 8 wells
K 7780	STD	Standards, ready-to-use (0, 0.1, 0.3, 0.6, 1.5, 4.0 μ M)	6 x 500 μ l
K 7780	CTRL 1	Control, ready-to-use (see specification for range)	1 x 500 μ l
K 7780	CTRL 2	Control, ready-to-use (see specification for range)	1 x 500 μ l
K 0006.C.100	WASHBUF A	Wash buffer concentrate, 10x	2 x 100 ml
K 7780	AB	SDMA antibody, ready-to-use	1 x 6 ml
K 7780	CONJ	Conjugate, ready-to-use	1 x 12 ml
K 0012.15	DERBUF	Reaction buffer, ready-to-use	1 x 15 ml
K 7780	DER	Derivatisation reagent, ready-to-use	1 x 6 ml
K 0002.15	SUB	Substrate (tetramethylbenzidine), ready-to-use	1 x 15 ml
K 0003.15	STOP	Stop solution, ready-to-use	1 x 15 ml

For reorders of single components, use the catalogue number followed by the label as product number.

4. MATERIAL REQUIRED BUT NOT SUPPLIED

- Ultra pure water*
- Calibrated precision pipets and 10-1000 μ l tips
- Foil to cover the microtiter plate
- Horizontal microtiter plate shaker
- Multi-channel pipets or repeater pipets
- Vortex
- Centrifuge, 3000 g
- Standard laboratory glass or plastic vials, cups, etc.
- Microtiter plate reader (required filters see chapter 7)

* Immundiagnostik AG recommends the use of Ultra Pure Water (Water Type 1; ISO 3696), which is free of undissolved and colloidal ions and organic molecules (free of particles > 0.2 μ m) with an electrical conductivity of 0.055 μ S/cm at 25 °C (\geq 18.2 M Ω cm).

5. STORAGE AND PREPARATION OF REAGENTS

- To run the assay more than once, ensure that reagents are stored at the conditions stated on the label.
- **Preparation of the wash buffer:** The **wash buffer concentrate (WASHBUF A)** has to be diluted with ultra pure water **1:10** before use (100 ml WASHBUF A + 900 ml ultra pure water), mix well. Crystals could occur due to high salt concentration in the concentrate. Before dilution, the crystals have to be redissolved at room temperature or in a water bath at 37 °C. The **WASHBUF A** is stable at **2-8 °C** until the expiry date stated on the label. **Wash buffer** (1:10 diluted WASHBUF A) can be stored in a closed flask at **2-8 °C for 1 month**.
- Store **standards and controls (STD/CTRL)** frozen at **-20 °C**. They are stable at -20 °C until the expiry date stated on the label. Thaw before use in the test and mix well. Re-freeze standards and controls after use.
- The ready-to-use **derivatisation reagent (DER)** is already dissolved in DMSO and is stable at **2-8 °C** until the expiry date stated on the label. DMSO crystallises at 2-8 °C. Before opening the DER, bring to room temperature and ensure that all crystals are dissolved. Please note: DMSO attacks all plastics but not polypropylene products and laboratory glass.
- All other test reagents are ready-to-use. Test reagents are stable until the expiry date (see label) when stored at **2-8 °C**.

6. STORAGE AND PREPARATION OF SAMPLES

Serum, EDTA and Li-heparin plasma

- Freshly collected serum, EDTA and Li-heparin plasma samples can be stored for up to 48 hours at room temperature or for 3 days at 2-8 °C. For longer storage keep samples frozen at -20 °C.
- The samples are analysed **undiluted**.
If the sample volume is less than 50 µl, we recommend a 1:2 dilution in reaction buffer (25 µl sample + 25 µl DERBUF). This dilution factor must be considered in data evaluation.
- For sample preparation, a derivatisation reagent for derivatisation of SDMA is added (see sample preparation procedure).

7. ASSAY PROCEDURE

Principle of the test

This ELISA is designed for the quantitative determination of SDMA. The assay is based on the method of competitive enzyme linked immunoassays.

The sample preparation includes the addition of a derivatisation reagent for SDMA derivatisation. Afterwards, the treated samples and the polyclonal SDMA antiserum are incubated in wells of a microtiter plate coated with SDMA derivative (tracer). During the incubation period, the target SDMA in the sample competes with the tracer, immobilised on the wall of the microtiter wells, for the binding of the polyclonal antibodies.

During the second incubation step, a peroxidase conjugated antibody is added to detect the anti-SDMA antibodies. After washing away the unbound components, tetramethylbenzidine (TMB) is added as a peroxidase substrate. Finally, the enzymatic reaction is terminated by an acidic stop solution. The colour changes from blue to yellow and the absorbance is measured in a photometer at 450 nm. The intensity of the yellow colour is inverse proportional to the SDMA concentration in the sample; this means high SDMA concentration in the sample reduces the concentration of tracer-bound antibodies and lowers the photometric signal. A dose response curve of absorbance unit (optical density, OD at 450 nm) vs. concentration is generated using the values obtained from the standards. SDMA, present in the patient samples, is determined directly from this curve.

Sample preparation procedure

Bring **all reagents and samples to room temperature** (15-30 °C) and mix well.

Derivatisation of standards, controls and samples is carried out in single analysis in vials (e.g. 1.5 ml vials).

We recommend preparing one derivatisation per standard, control and sample and transferring it in duplicate determinations into the wells of the microtiter plate.

1.	Add 200 µl standard (STD), 200 µl control (CTRL) and 50 µl sample in the corresponding vials.
2.	Add 150 µl reaction buffer (DERBUF) only to the samples .

3.	Add 50 µl derivatisation reagent into each vial (STD, CTRL, sample), mix thoroughly by repeated inversion or several seconds on a vortex mixer. Incubate for 45 min at room temperature (15-30 °C) on a horizontal shaker .
----	---

2 x 50 µl of the derivatised standards, controls and samples are used in the ELISA as duplicates.

Test procedure

Mark the positions of standards/controls/samples in duplicate on a protocol sheet.

Take as many microtiter strips as needed from the kit. Store unused strips covered at 2-8 °C. Strips are stable until expiry date stated on the label.

4.	For the analysis in duplicate, take 2 x 50 µl of the derivatised standards/ controls/samples out of the vials and add into the respective wells of the microtiter plate.
5.	Add 50 µl SDMA antibody (AB) into each well of the microtiter plate.
6.	Cover the strips and incubate for 2 hours at room temperature (15-30 °C) on a horizontal shaker .
7.	Discard the content of each well and wash 5 times with 250 µl wash buffer . After the final washing step, remove residual wash buffer by firmly tapping the plate on absorbent paper.
8.	Add 100 µl conjugate (CONJ) into each well.
9.	Cover the strips and incubate for 1 hour at room temperature (15-30 °C) on a horizontal shaker .
10.	Discard the content of each well and wash 5 times with 250 µl wash buffer . After the final washing step, remove residual wash buffer by firmly tapping the plate on absorbent paper.
11.	Add 100 µl substrate (SUB) into each well.
12.	Incubate for 10-15 min* at room temperature (15-30 °C) in the dark .
13.	Add 100 µl stop solution (STOP) into each well and mix well.

- | | |
|-----|--|
| 14. | Determine absorption immediately with an ELISA reader at 450 nm against 620 nm (or 690 nm) as a reference. If no reference wavelength is available, read only at 450 nm. If the extinction of the highest standard exceeds the range of the photometer, absorption must be measured immediately at 405 nm against 620 nm (690 nm) as a reference. |
|-----|--|

* The intensity of the colour change is temperature sensitive. We recommend observing the colour change and stopping the reaction upon good differentiation.

For automated ELISA processors, the given protocol may need to be adjusted according to the specific features of the respective automated platform. For further details please contact your supplier or Immundiagnostik AG.

8. RESULTS

The following algorithms can be used alternatively to calculate the results. We recommend using the 4 parameter algorithm.

1. 4 parameter algorithm

It is recommended to use a linear ordinate for optical density and a logarithmic abscissa for concentration. When using a logarithmic abscissa, the zero standard must be specified with a value less than 1 (e.g. 0.001).

2. Point-to-point calculation

We recommend a linear ordinate for optical density and a linear abscissa for concentration.

3. Spline algorithm

We recommend a linear ordinate for optical density and a linear abscissa for concentration.

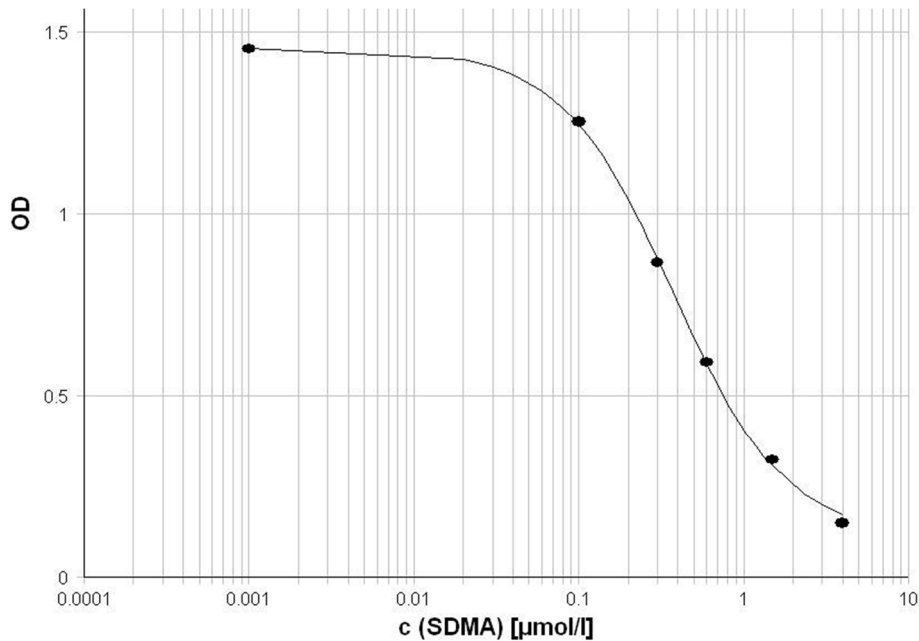
The plausibility of the pairs of values should be examined before the automatic evaluation of the results. If this option is not available with the used program, a control of the paired values should be done manually.

Serum, EDTA and Li-heparin plasma

The concentrations can be determined directly from the standard curve in $\mu\text{mol/l}$. **No factor** is required.

Consider the dilution factor if samples have been pre-diluted 1:2.

In the following, an example of a calibration curve is given; do not use it for the calculation of your results.



9. LIMITATIONS

Samples with an OD lower than the OD of the highest standard should be diluted with reaction buffer and re-assayed. Please consider this dilution factor when calculating the results.

10. QUALITY CONTROL

Immundiagnostik recommends the use of external controls for internal quality control, if possible.

Control samples should be analysed with each run. Results, generated from the analysis of control samples, should be evaluated for acceptability using appropriate statistical methods. The results for the patient samples may not be valid if within the same assay one or more values of the quality control samples are outside of the acceptable limits.

Reference range

Based on internal studies with human serum samples from apparently healthy persons ($n = 40$), a mean value of $0.47 \mu\text{mol/l}$ was estimated. The standard deviation (SD) was $0.09 \mu\text{mol/l}$. From mean value ± 2 SD a normal range of $0.29 - 0.65 \mu\text{mol/l}$ was estimated.

We recommend each laboratory to establish its own reference range.

11. PERFORMANCE CHARACTERISTICS

Precision and reproducibility

Intra-assay (n = 10)

Sample	SDMA [$\mu\text{mol/l}$]	CV [%]
1	0.27	7.5
2	0.67	4.8

Inter-assay (n = 6)

Sample	SDMA [$\mu\text{mol/l}$]	CV [%]
1	0.22	9.9
2	0.63	7.4

Spiking recovery

Human

One serum sample was spiked with different SDMA concentrations and measured in this assay. The mean recovery rate was 100.4 % (n = 5).

Spike [$\mu\text{mol/l}$]	SDMA expected [$\mu\text{mol/l}$]	SDMA measured [$\mu\text{mol/l}$]	Recovery [%]
0		0.736	
0.5	1.236	1.220	98.7
1	1.736	1.774	102.2

Dog

Two serum samples were spiked with different SDMA concentrations. The mean recovery rate was 94.1 % (n = 2).

Spike [$\mu\text{mol/l}$]	SDMA expected [$\mu\text{mol/l}$]	SDMA measured [$\mu\text{mol/l}$]	Recovery [%]
0		0.526	
0.5	1.026	0.937	91.3
1	1.526	1.561	102.3
0		0.454	
0.5	0.954	0.897	94.0
1	1.454	1.288	88.6

Cat

Two serum samples were spiked with different SDMA concentrations. The mean recovery rate was 95.4 % (n = 2).

Spike [μmol/l]	SDMA expected [μmol/l]	SDMA measured [μmol/l]	Recovery [%]
0		0.395	
0.5	0.895	0.833	93.1
1	1.395	1.252	89.8
0		0.326	
0.5	0.826	0.818	99.0
1	1.326	1.321	99.6

*Dilution recovery***Human**

One spiked sample was diluted with reaction buffer. The mean recovery rate was 90.4 % (n = 5).

Dilution	SDMA expected [μmol/l]	SDMA measured [μmol/l]	Recovery [%]
		1.774	
1:2	0.887	0.876	98.8
1:4	0.444	0.364	82.1

Dog

Two samples were diluted with serum containing 0.126 μmol/l SDMA. The mean recovery rate was 109.0 % (n = 2).

Dilution	SDMA expected [μmol/l]	SDMA measured [μmol/l]	Recovery [%]
		0.526	
1:2	0.326	0.334	102.5
1:4	0.226	0.241	106.6
		0.454	
1:2	0.290	0.340	117.2
1:4	0.208	0.228	109.6

Cat

Two samples were diluted with serum containing 0.126 µmol/l SDMA. The mean recovery rate was 97.1 % (n = 2).

Dilution	SDMA expected [µmol/l]	SDMA measured [µmol/l]	Recovery [%]
		0.395	
1:2	0.261	0.247	94.8
1:4	0.193	0.194	100.4
		0.326	
1:2	0.226	0.231	102.2
1:4	0.176	0.160	90,9

Analytical sensitivity

The zero-standard (STD 1) was measured 40 times. The detection limit was set as $B_0 - 2 \text{ SD}$ and estimated to be 0.05 µmol/l.

Sample	Mean value [OD]	2 x standard deviation (2 SD)	Detection limit [µmol/l]
zero-standard	2.3	0.14	0.05

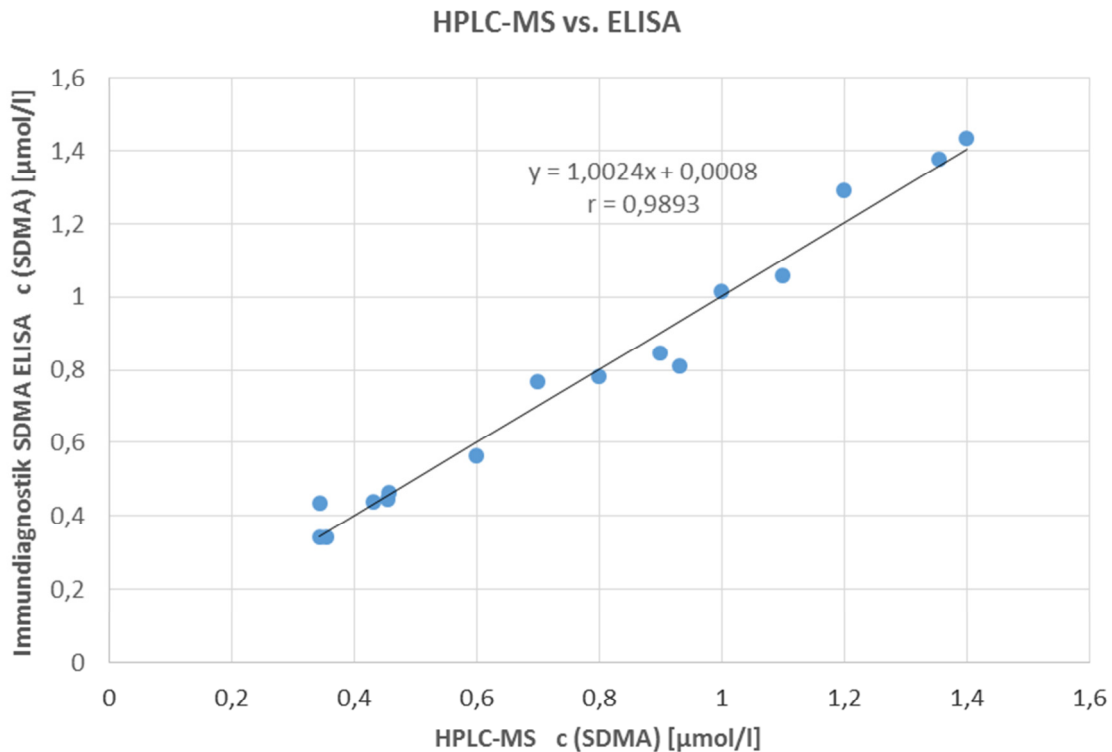
Specificity

The specificity of the antibody was tested by measuring the cross-reactivity against a range of compounds with structural similarity to SDMA. The specificity is calculated in percent in relation to the SDMA-binding activity.

ADMA	< 0.1 %
L-arginine	< 0.001 %

Correlation with HPLC-MS

16 samples were measured with this ELISA and HPLC-MS. The correlation was $r = 0.99$.



12. PRECAUTIONS

- All reagents in the kit package are for *in vitro* diagnostic use only.
- Human materials used in kit components were tested and found to be negative for HIV, Hepatitis B and Hepatitis C. However, for safety reasons, all kit components should be treated as potentially infectious.
- Kit reagents contain sodium azide or ProClin as bactericides. Sodium azide and ProClin are toxic. Substrates for the enzymatic colour reactions are toxic and carcinogenic. Avoid contact with skin or mucous membranes.
- The stop solution consists of diluted sulfuric acid, a strong acid. Although diluted, it still must be handled with care. It can cause burns and should be handled with gloves, eye protection, and appropriate protective clothing. Any spill should be wiped up immediately with copious quantities of water. Do not breathe vapour and avoid inhalation.

13. TECHNICAL HINTS

- Do not interchange different lot numbers of any kit component within the same assay. Furthermore, we recommend not assembling wells of different microtiter plates for analysis, even if they are of the same batch.
- Control samples should be analysed with each run.
- Reagents should not be used beyond the expiration date stated on the kit label.
- Substrate solution should remain colourless until use.
- To ensure accurate results, proper adhesion of plate sealers during incubation steps is necessary.
- Avoid foaming when mixing reagents.
- Do not mix plugs and caps from different reagents.
- The assay should always be performed according to the enclosed manual.

14. GENERAL NOTES ON THE TEST AND TEST PROCEDURE

- This assay was produced and distributed according to the IVD guidelines of 98/79/EC.
- The guidelines for medical laboratories should be followed.
- Incubation time, incubation temperature, and pipetting volumes of the components are defined by the producer. Any variation of the test procedure, which is not coordinated with the producer, may influence the results of the test. Immundiagnostik AG can therefore not be held responsible for any damage resulting from incorrect use.
- Warranty claims and complaints regarding deficiencies must be logged within 14 days after receipt of the product. The product should be sent to Immundiagnostik AG along with a written complaint.

15. REFERENCES

General literature












1. Fleck C, Schweitzer F, Karge E, Busch M, Stein G. Serum concentrations of asymmetric (ADMA) and symmetric (SDMA) dimethylarginine in patients with chronic kidney diseases. *Clinical Chimica Acta* **336**: 1 – 12 (2003).

2. D'Apolito O, Paglia G, Tricarico F, Garofalo D, Pilotti A, Lamacchia O, Cignarelli M, Corso G. Development and validation of a fast quantitative method for plasma dimethylarginines analysis using liquid chromatography-tandem mass spectrometry. *Clinical Biochemistry* **41**: 1391 – 1395 (2008).
3. Kielstein JT, Salpeter SR, Bode-Böger SM, Cooke JP, Fliser D. Symmetric dimethylarginine (SDMA) as endogenous marker of renal function – a meta-analysis. *Nephrol. Dial. Transplant* **21**: 2446 - 2451(2006).
4. Bode-Böger SM, Scalera F, Kielstein JT, Martens-Lobenhoffer J, Breithardt G, Fobker M, Reinecke H. Symmetrical Dimethylarginine: A new combined parameter for renal function and extent of coronary artery disease. *J. Am. Soc. Nephrol.* **17**: 1128 – 1134 (2006).

Literature using Immundiagnostik SDMA ELISA

5. Koch A, Weiskirchen R, Bruensing J, Dückers H, Buendgens L, Kunze J, Matthes M, Luedde T, Trautwein C, Tacke F. Regulation and prognostic relevance of symmetric dimethylarginine serum concentrations in critical illness and sepsis. *Mediators of Inflammation* **2013**: 413826 (2013).
6. Tenderenda-Banasiuk E, Wasilewska A, Taranta-Janusz K, Korzeniecka-Kozerska A. Asymmetric and symmetric dimethylarginine in adolescents with hyperuricemia. *Disease Markers* **35**: 407-412 (2013).

Used symbols:

	Temperature limitation		Catalogue Number
	In Vitro Diagnostic Medical Device		To be used with
	Manufacturer		Contains sufficient for <n> tests
	Lot number		Use by
	Attention		Consult instructions for use
	Consult specification data sheet		