



Ultra-Sensitive Measurement of Methotrexate

Methotrexate ELISA Kit

A quantitative approach to monitoring methotrexate levels during treatment of cancer and autoimmune disease

Methotrexate ELISA Kit (ENZ-KIT142)

Methotrexate is a drug used in the treatment of cancer and autoimmune disease. It is designed as an anti-folate to inhibit the metabolism of folic acid. Monitoring methotrexate levels is important to assure appropriate levels are maintained during therapy or treatment. High levels of methotrexate can lead to toxicity and potential renal failure as well as immunosuppression. Additionally, methotrexate is known to interact with a wide variety of drugs leading to additional complications. Determining the presence of methotrexate in samples from subjects in blinded research studies can assist in the interpretation of study results.

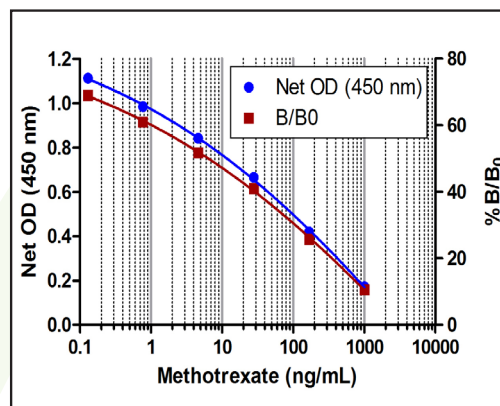
Key Features

- Sensitive measurement of methotrexate, detecting as little as 0.087 ng/ml
- Broad dynamic range minimizing the need for dilution
- Compatible with various samples including serum, plasma and urine in human, mouse, and rat samples
- High throughput format with results in 1.5 hours for up to 38 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Research Areas

*Cancer
Rheumatoid Arthritis
Inflammatory Bowel Disease
Psoriasis
Bioprocess*

Ultra-Sensitive Measurement of Methotrexate



Standard curve experiment using Methotrexate ELISA Kit (ENZ-KIT142)

PRODUCT NAME	PRODUCT #
Methotrexate ELISA Kit	ENZ-KIT142
Methotrexate, pAb	ADI-905-052
Methotrexate	ALX-440-045

Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com