



Sensitive Measurement of MBL

MBL Oligomer ELISA Kit

High throughput format with results in just 4 hours

MBL Oligomer ELISA Kit (BDP-KIT-029)

MBL (Mannan-Binding Lectin) is a carbohydrate-binding protein produced in the liver and secreted into the blood. The protein plays a key factor in innate immunity, functioning as a player in the complement activation cascade and defending the body against invading microorganisms (bacteria, viruses, protozoa and fungi). The MBL Oligomer ELISA kit is for the measurement of oligomerized MBL in human samples. Absorbance is read at 450nm. Mannan-binding lectin (MBL) is an important component of the innate immune system.

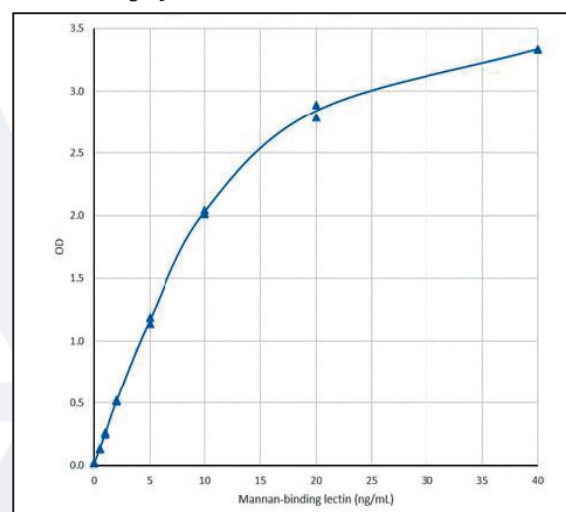
Benefits

- Highly sensitive measurement, detecting as low of 2 ng/ml of oligomerized MBL
- High throughput format with results in just 4 hours for up to 40 samples in duplicate
- Ready-to-use pre-coated plate and individual calibrators saves time and reduce errors
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Research Areas

- Cancer
- Cardiovascular
- Dengue
- Diabetes
- Hepatitis B
- HIV
- Immunology
- Pulmonary Fibrosis
- Renal Disease
- Rheumatoid Arthritis

Highly-sensitive Measurement of MBL



Typical standard curve for MBL Oligomer ELISA Kit (BDP-KIT-029)

RELATED PRODUCTS	PRODUCT #
CD40L, Soluble (human) ELISA Kit (high sensitivity)	ALX-850-311
Complement C3a des Arg (human) ELISA Kit	ADI-900-058
Complement C4a des Arg (human) ELISA Kit	ADI-900-059
COX-2 (human) ELISA Kit	ADI-900-094
CRP ELISA Kit	ENZ-KIT102
FasL, Soluble (human) ELISA Kit	ALX-850-246
MBL (human) mAb	BPD-HYB-131
MEGACD40L® Protein (soluble)(human)(recombinant)	ALX-522-110
NFκB p65 ELISA Kit (chemiluminescent)	ADI-EKS-446
SUPERFAS LIGAND® (soluble)(human)(recombinant)	ALX-522-020

Fuel 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