



Highly Sensitive ELISA for Inflammation Research

CRP ELISA Kit

Highly sensitive ELISA kit for the quantification of CRP, an early biomarker for inflammation

CRP ELISA Kit (ENZ-KIT102)

C-Reactive Protein (CRP) is a pentraxin family protein that is widely used as a biomarker for systemic inflammation and tissue injury. Its primary role in innate immunity is to function as an opsonin for pathogens via activation of the complement pathway and through binding to Fc gamma receptors. CRP has been shown to play an important role in a variety of pathological pathways such as chronic inflammatory disorders, myocardial infarction, pulmonary disorders, atherosclerosis, cancer, metabolic syndrome, and depression.

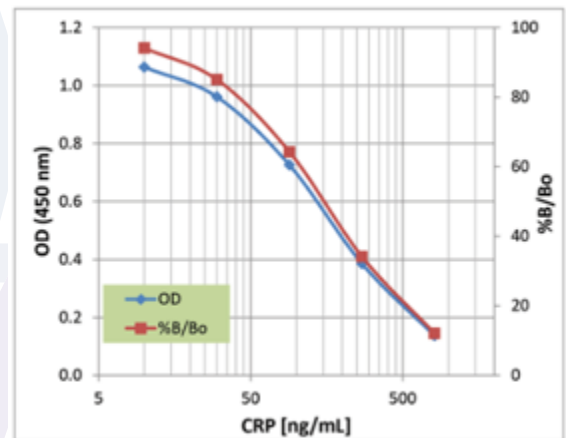
Key Features

- Sensitive measurement of CRP, detecting as little as 12.685 ng/ml
- Negligible reactivity for similar family members
- Compatible with plasma and serum human samples
- High throughput format with results in < 2 hours for up to 40 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Research Areas

Cancer *Inflammation*
Cardiovascular *Metabolic Syndrome*
Depression *Pulmonary*

High Sensitivity Measurement of CRP



Standard curve experiment using CRP ELISA Kit (ENZ-KIT102)

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
CRP (human), ELISA Kit	ENZ-KIT102	IL-1F8 (human), (recombinant)	ALX-522-129-C050
Complement C3a des Arg (human), ELISA Kit	ADI-900-058	IL-1F9 (human), (recombinant)	ALX-522-130-C050
Complement C4a des Arg (human), ELISA Kit	ADI-900-059	IL-6 (human), ELISA Kit	ADI-900-033
COX-2 (human), ELISA Kit	ADI-900-094	IL-8 (human), ELISA Kit	ADI-900-156
C-Reactive Protein (human), (recombinant)	ENZ-PRT164	IL-33 (human), ELISA Kit	ADI-900-201
Haptoglobin (human), ELISA Kit	ADI-900-229-0001	PGE2 ELISA Kit	ADI-930-001
Histamine ELISA Kit	ENZ-KIT140	PGE2 High Sensitivity ELISA Kit	ADI-930-001
IL-1β (soluble) (human), (recombinant)	ALX-522-056-C010	TNF-α (human), ELISA Kit	ADI-900-099
IL-1β (human), ELISA Kit	ADI-900-130	TNF-α (soluble) (human), (recombinant)	ALX-522-008-C050
IL-1F6 (human), (recombinant)	ALX-522-128-C050	Trefoil Factor 1 (human), ELISA Kit	ALX-850-382-KI01

Activate 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