



SMN Detection in 3 Hours

SMN ELISA Kit (ADI-900-209)

First-to-market SMN protein immunoassay

Survival Motor Neuron (SMN) is a ~38 kDa protein produced chiefly by the SMN1 gene. A nearly identical centromeric copy of the gene (SMN2) also produces a small amount of full-length SMN protein, but due to a translationally silent C-T transition that results in alternative splicing of the pre-mRNA, most of the resulting SMN is truncated, causing reduced protein stability and lower overall SMN levels. Deletion or mutation of the SMN1 gene results in a reduced level of full-length SMN protein and manifests as a range of neuromuscular phenotypes in humans as the disease spinal muscular atrophy (SMA). SMA is characterized by muscle weakness and atrophy, functional disability and is the most common lethal genetic disease of infants and toddlers. The SMN ELISA Kit was developed in collaboration with the SMA Foundation in New York City.

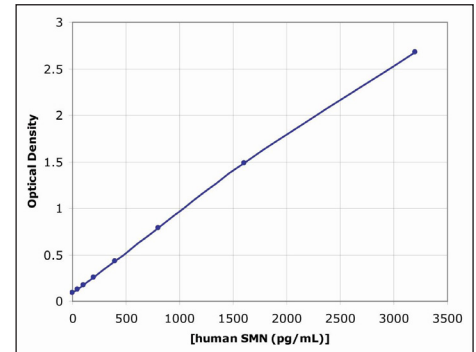
Key Features

- Highly sensitive, measure as little as 50 pg/mL of SMN
- Compatible with human and mouse SMN in cell lysate samples
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Results from up to 39 samples in duplicate in just <3 hours
- Ready-to-use liquid color coded reagents and pre-coated 96-well plate

Areas of Research

- Aging
- RNA Metabolism
- Neuroscience
- SMA

Highly Sensitive Measurement of SMN



The SMN ELISA kit is a colorimetric immunometric immunoassay kit with results in 3 hours.

RELATED PRODUCTS

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
HSP25, pAb	ADI-SPA-801	IMMUNOSETS® HSP25 (rodent), ELISA Development Set	ADI-960-075
HSP27 (human), ELISA Kit	ADI-EKS-500	[pSer ¹⁵]HSP27, pAb	ADI-SPA-525
HSP27 (human), (rec.)	ADI-SPP-715	[pSer ⁷⁸]HSP27, pAb	ADI-SPA-523
HSP27, mAb (G3.1)	ADI-SPA-800	[pSer ⁸²]HSP27, pAb	ADI-SPA-524
HSP27, pAb	ADI-SPA-803	SMN1 (human), (rec.) (His-tag)	ADI-NBP-201
IMMUNOSETS® HSP27 high sensitivity (human), ELISA development set	ADI-960-076	SMN1, mAb (2B1)	ADI-NBA-201

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