

SMN ELISA kit

First ready-to-use SMN ELISA kit commercially available for neurodegenerative disease research.

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

Survival Motor Neuron (SMN) is a ~38 kDa protein produced chiefly by the SMN1 gene, located on the telomeric portion of chromosome 5q. 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. Approximately one in 35 adults is a carrier of the SMN1 mutation. The incidence of SMA is 1 in 6,000 to 1 in 10,000 live births. SMN protein is present in the cell cytoplasm, and also in the nucleus where it is concentrated in "gem" structures associated with Cajal bodies. SMN protein is a constituent of Gemin-containing complexes, and is thought to participate in many aspects of RNA metabolism. SMN complexes have been shown to mediate the assembly of uridine rich small nuclear ribonucleoproteins (snRNPs), which in turn act as critical components of spliceosomes.

Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

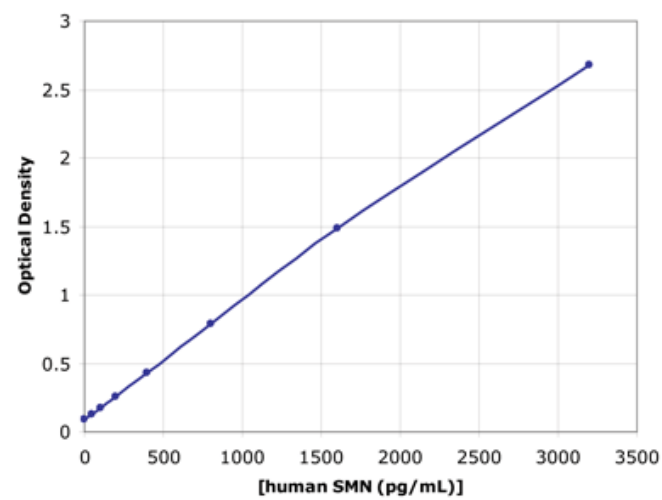
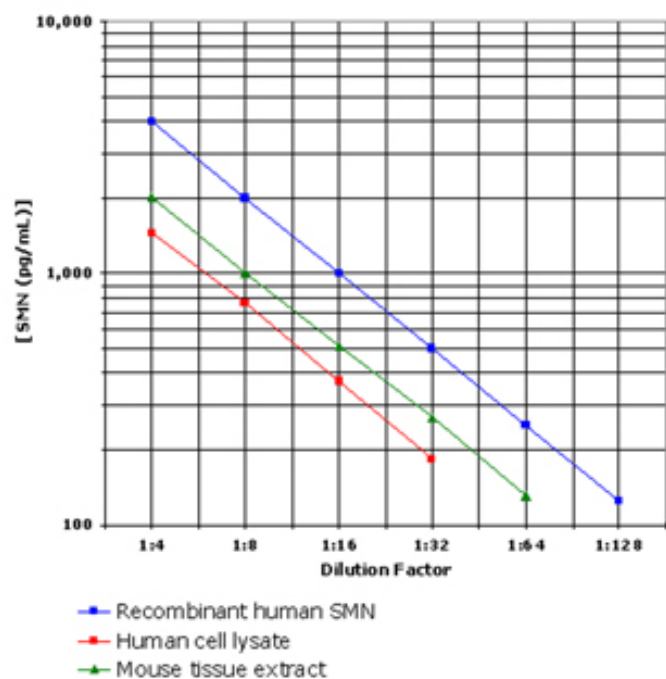
ADI-900-209

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement, detecting as little as 50 pg/mL of SMN
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Convenient and ready-to-use liquid color coded reagents and pre-coated 96-well plate to save time and minimize errors
- High throughput format with results in 3 hours for up to 39 samples in duplicate



Handling & Storage

Use/Stability Store all components at +4°C, except Standard at -20°C.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Survival motor neuron

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human and mouse SMN in cell lysate samples.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter Plate, Conjugate, Antibody, Assay Buffer 13, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Extraction Reagent 4

Crossreactivity No cross reactivity.

Sensitivity 50pg/ml (range 50-3200pg/ml)

Species Reactivity Human, Mouse

Technical Info / Product Notes The kit was developed in collaboration with the SMA Foundation (New York, NY).
Please read the complete kit insert before performing this assay.
Protected by US Patent no. US. 6,080,577.

UniProt ID Q16637

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)