

24(S)- Hydroxycholesterol ELISA kit

Only quantitative 24(S)-hydroxycholesterol ELISA kit available for neurodegenerative research

The 24(S)-Hydroxycholesterol ELISA kit is a complete, colorimetric, competitive immunoassay kit for the quantitative determination of 24(S)-Hydroxycholesterol in tissue culture media (culture supernatant), cerebral spinal fluid and tissue homogenate samples with results in just 2 hours.

The homeostasis and trafficking of cholesterol is an essential component of both the central and peripheral nervous system in the maintenance of neuronal tissues, and disturbances in this homeostasis may be due to the onset of various neurological diseases such as Alzheimer's Disease, Huntington's Disease and multiple sclerosis. Apolipoprotein E and Cyp46 (also known as 24S-Cholesterol Hydroxylase) are both important in the homeostasis of cerebral cholesterol and thus are of clinical interest in understanding the relation of these molecules with the pathogenesis of these neurodegenerative diseases.

24-OHC, an enzymatically-generated side chain-hydroxylated derivative of cholesterol, is a pivotal marker in the study of cerebral cholesterol homeostasis. Cholesterol is unable to cross the blood-brain barrier however, Cyp46 enzyme converts cholesterol to the more soluble 24-OHC, and this hydroxylated form of cholesterol is able to cross the blood-brain barrier. This conversion allows for the reduction of cholesterol in the brain and the efflux of 24-OHC from the brain into cerebral spinal fluid and blood. The flux of 24-OHC has been seen in patients with a variety of neurodegenerative diseases. In the instance of Alzheimer's disease, the change in 24S-hydroxycholesterol concentrations may be indicative of different pathogenetic mechanisms and/or the progression of the disease. As in the case of multiple sclerosis, concentrations of 24-OHC have been shown to decrease, likely due to the loss of neuronal cells responsible for the synthesis.

Citations: 11

[View Online »](#)

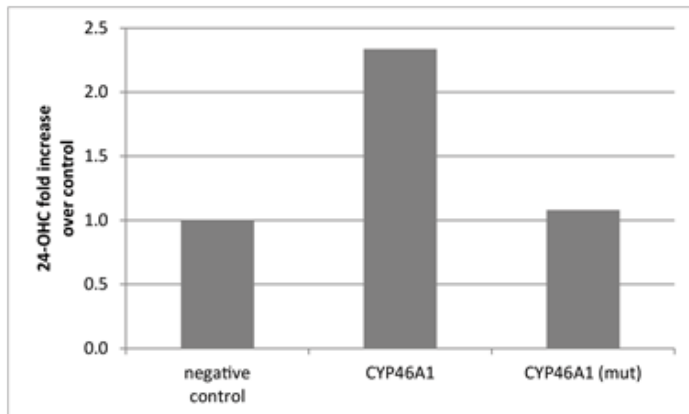
Ordering Information

[Order Online »](#)

ADI-900-210-0001

96 wells

- Convenient – user-friendly alternative to mass spectrometry
- Sensitive – measure as little as 0.78 ng/ml of 24(S)-hydroxycholesterol
- Time Saving – results from up to 36 samples in duplicate in just 2 hours
- Specific – low cross-reactivity with structurally related molecules
- Easy-to-use – liquid color-coded reagents reduce errors



Download our Application Note: [A Competitive, Colorimetric ELISA for Quantification of 24\(S\)-Hydroxycholesterol, a Biomarker for Neurodegeneration](#)

24(S)-Hydroxycholesterol ELISA kit 24-OHC-fold-increase, 24(S)-Hydroxycholesterol ELISA kit 24-OHC-fold-increase, yes

ADI-900-210, ADI-900-210-

parallelism.gif, <http://static.enzolifesciences.com/fileadmin/files/image/ADI-900-210-parallelism.gif>, 24(S)-Hydroxycholesterol

ELISA kit Parallelism curve, 24(S)-Hydroxycholesterol

ELISA kit Parallelism curve, yes

ADI-900-210, ADI-900-

210_std_curve.jpg, http://static.enzolifesciences.com/fileadmin/files/image/ADI-900-210_std_curve.jpg, 24(S)-Hydroxycholesterol

ELISA kit Standard curve, 24(S)-Hydroxycholesterol

ELISA kit Standard curve, yes

ADI-900-210, ADI-900-210-

experiment_eval.gif, http://static.enzolifesciences.com/fileadmin/files/image/ADI-900-210-experiment_eval.gif, Detect normal levels of

24(S)-hydroxycholesterol. Normal human cerebral spinal fluid samples were diluted 1:2 in assay buffer and analyzed in the assay for 24(S)-Hydroxycholesterol levels. The blue bars represent the levels of 24(S)-Hydroxycholesterol in each sample as determined by ELISA. The dashed lines represent the maximum and minimum of the range of endogenous CSF 24(S)-Hydroxycholesterol levels as reported in the literature (Annu. Rev. Biochem. 2009.78:1017-1040.), 24(S)-Hydroxycholesterol ELISA kit experiment_eval, 24(S)-Hydroxycholesterol ELISA kit experiment_eval, yes

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, http://static.enzolifesciences.com/fileadmin/files/image/ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

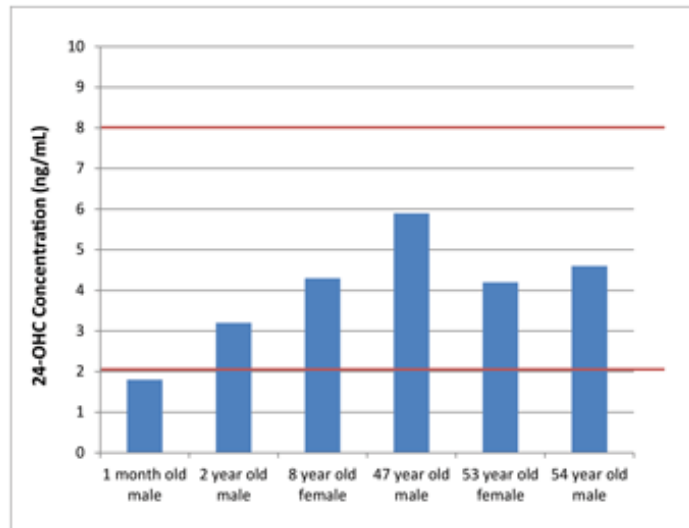
ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

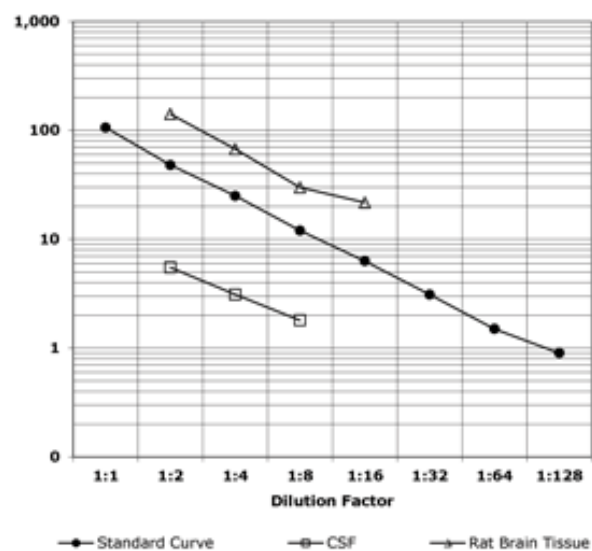
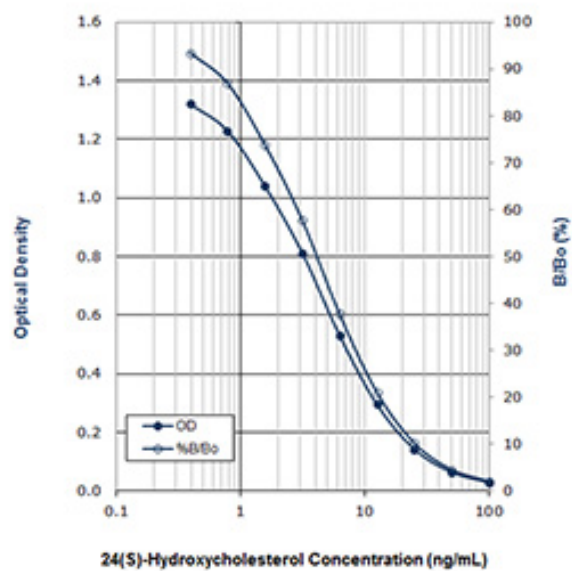
ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)

ADI-900-212, ADI-900-212_ELISA-WB_01.jpg, Correlation of p62 (Prod. No. ADI-900-212) and NBR1 (Prod. No. ADI-900-211) immunoassays to autophagy induction. MDA-MB-231 human breast cancer cells were treated with 2µM of withaferin A (WA)



Detect normal levels of 24(S)-hydroxycholesterol. Normal human cerebral spinal fluid samples were diluted 1:2 in assay buffer and analyzed in the assay for 24(S)-Hydroxycholesterol levels. The blue bars represent the levels of 24(S)-Hydroxycholesterol in each sample as determined by ELISA. The dashed lines represent the maximum and minimum of the range of endogenous CSF 24(S)-Hydroxycholesterol levels as reported in the literature (Annu. Rev. Biochem. 2009.78:1017-1040.)



Handling & Storage

Use/Stability All kit components should be stored at 4°C.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of species independent 24(S)-Hydroxycholesterol in cerebral spinal fluid, culture supernatants and tissue samples.

Assay Time 2 hours

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

Contents Microtiter Plate, Conjugate, Antibody, Assay Buffer, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Streptavidin-HRP

Sensitivity 0.78 ng/ml (range 0.39-100ng/ml)

Species Reactivity Species independent

Wavelength 450 nm

Last modified: May 29, 2024



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)