

LVV Hemorphin 7

ELISA kit

Most sensitive and rapid LVV Hemorphin 7 kit for neuroscience market.

The LVV Hemorphin 7 EIA kit is a colorimetric, competitive immunoassay kit with results in just 3 hours.

Hemorphins are opioid peptides derived by proteolysis from hemoglobin. Their sequences are identical in several mammalian species including human, sheep and bovine. LVV hemorphin 7 (LVVYPWTQRF) binds strongly to the Angiotensin IV (AT4) receptors in the brain. The AT4 receptor is an integral membrane aminopeptidase also known as IRAP (insulin regulated membrane aminopeptidase). LVV hemorphin 7 and AT4 are not substrates but rather inhibitors of the AT4 (IRAP) receptor. Both promote learning and memory and reverse amnesia in animal models. Elevated serum levels of LVV Hemorphin 7 have also been documented in patients with some forms of breast cancers that are associated with an increased expression of cathepsins B and D.

Citations: 1

[View Online »](#)

Ordering Information

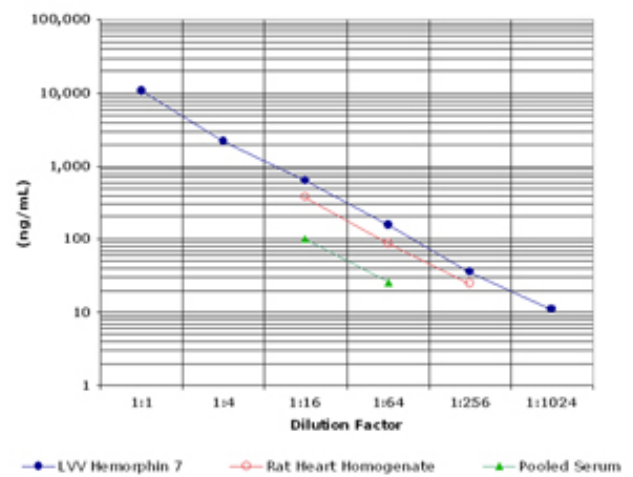
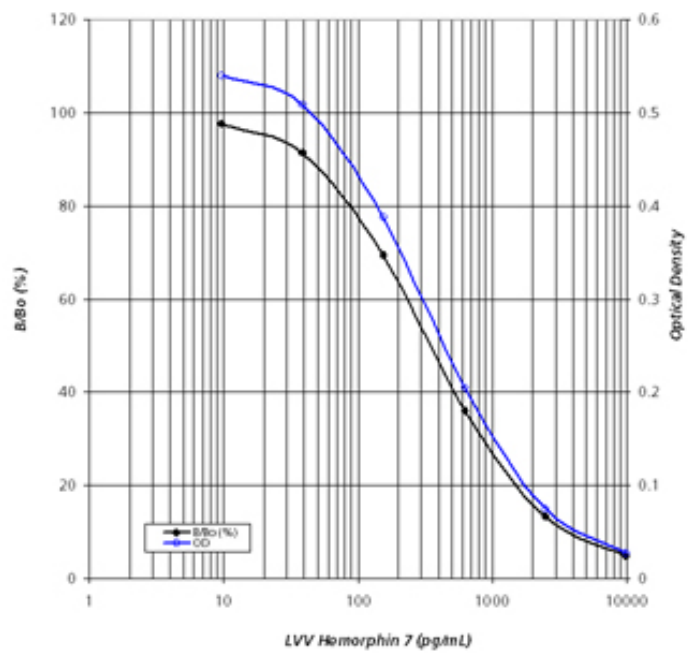
[Order Online »](#)

ADI-900-205	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - reliably measure as little as 6.1pg/ml of LVV Hemorphin 7
- Rapid - results in just 3 hours from up to 40 samples in duplicate
- Specific - low cross-reactivity to similar peptides
- Easy-to-use - liquid color-coded reagents reduce errors



Handling & Storage

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

Long Term Storage -20°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 LVV Hemorphin 7 in serum and tissue homogenate samples.

Assay Time 3 hours

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

Contents Microtiter Plate, Conjugate, Antibody, Assay Buffer 16, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Streptavidin-HRP

Crossreactivity LVV Hemorphin 7 (100%) and <0.003%: Leu-Valorphin-Arg (LVV Hemorphin-6), Valorphin, Angiotensin (1-12), Angiotensin I, Angiotensin (1-9), Angiotensin II, Angiotensin (1-7), Angiotensin A, Angiotensin III, Angiotensin IV

Sensitivity 6.1 pg/ml (range 9.8 – 10,000 pg/ml)

Species Reactivity Species independent

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)