

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 »](#)

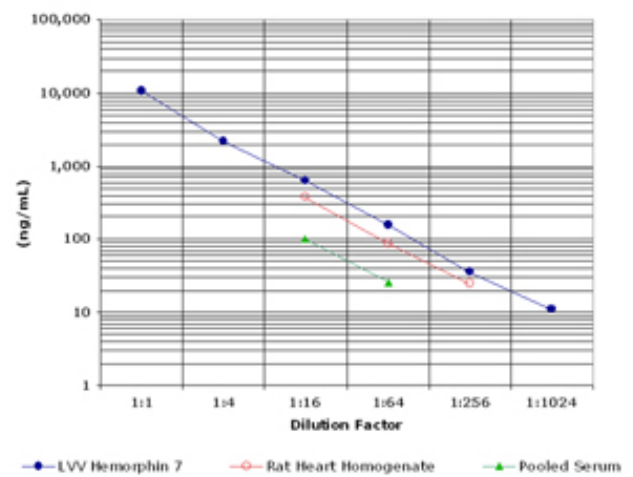
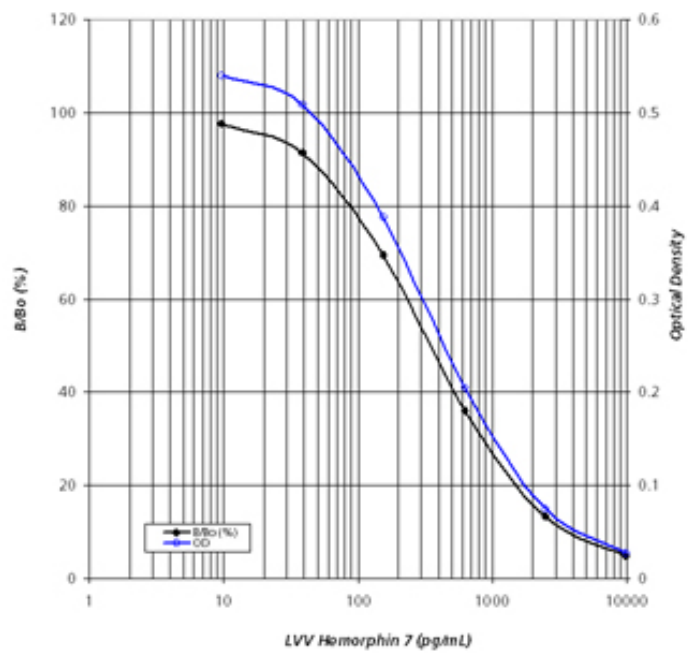
ADI-900-205

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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