

# Neurotensin 1 receptor (NTS<sub>1</sub>) polyclonal antibody

The subtype 1 neurotensin receptor (NTS1) is a seven-transmembrane domain containing G protein coupled receptor (GPCR) that mediates most of the known effects of neurotensin. NTS1 activates a phosphatidylinositol-calcium second messenger system in the central and peripheral nervous systems, mediating such functions as hypotension, hyperglycemia, hypothermia, antinociception, and the regulation of intestinal motility and secretion. NTS1 signaling is primarily mediated via coupling to Gq/G11 family G proteins, although coupling of NTS1 through Gs and Gi/o family proteins has been shown to modulate both adenylate and guanylate cyclase activity.

This antibody is covered by our [Worry-Free Guarantee](#).

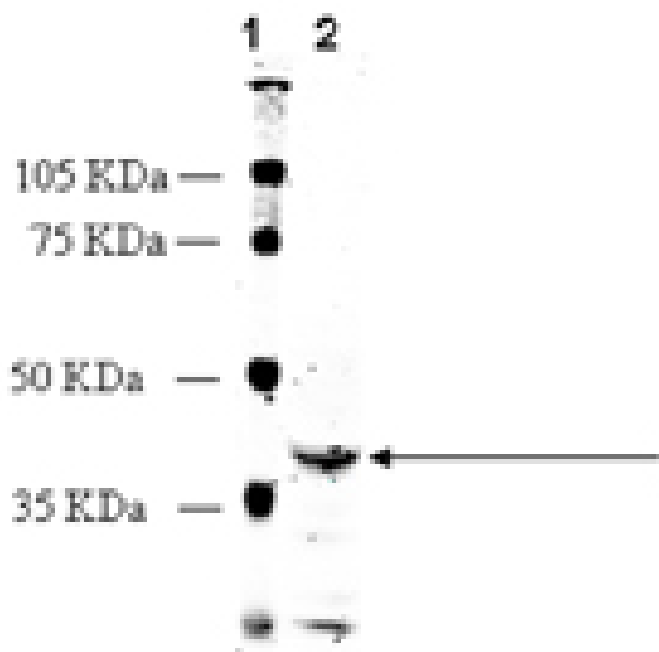
## Ordering Information

[Order Online »](#)

ADI-905-898-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW marker (1) and 20µg rat kidney extract (2) probed with Neurotensin Receptor 1 (NTS1) pAb.

## Handling & Storage

**Handling** Keep on ice at all times.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** IHC, WB

**Application Notes** Detects a band of ~40kDa. by Western blot.

**Crossreactivity** Predicted species reactivity based on epitope sequence homology: bovine, dog, rhesus macaque.

**Formulation** Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the N-terminus of human Neurotensin 1 Receptor (NTS1).

**Purity Detail** Peptide affinity purified.

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** P30989

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).

**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-be@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)