

Neurokinin-1 receptor (rat) polyclonal antibody

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

Ordering Information

[Order Online »](#)

BML-NA4300-0025	25µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NK1r, Substance P receptor
Application	IHC
Formulation	Liquid. Antiserum containing 0.1% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 393-407 (K393TMTTESSSFYSNMLA407) of C-terminal rat NK1r (neurokinin-1 receptor).
Source	From rabbit serum.
Species Reactivity	Rat
UniProt ID	P14600
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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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