FMRFamide polyclonal antibody

Note: Users should note that the antiserum is specific to the X-R-F-NH2 terminus and, accordingly, bioactive peptides with similar C-terminal sequences (e.g. mammalian g1-MSH9 and other N-terminally extended forms) may give strong positive immunostaining on certain tissues. The use of proteolytic enzymes (e.g. trypsin) to improve immunostaining should also be avoided as this may result in the appearance of 'false positive' staining due to the enzymatic induction of reactive -R-F-NH2 from nonreactive -R-F-G-K/R-K/R- pro-hormone sequences.

This antibody is covered by our Worry-Free Guarantee.

Citations: 6

View Online »

Ordering Information

Order Online »

BML-FA1155-0025	25µl
BML-FA1155-0100	100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Phe-Met-Arg-Phe-amide

Application IHC

Formulation Liquid. Neat serum containing 0.1% sodium azide.

Host Rabbit

Immunogen Synthetic FMRF-amide (Phe-Met-Arg-Phe-amide).

Source Whole rabbit serum.

Species Reactivity Species independent

Specificity Recognizes FMRF in multiple species including human.

UniProt ID O15130

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

