

# FMRamide

## polyclonal antibody

Note: Users should note that the antiserum is specific to the X-R-F-NH<sub>2</sub> 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-NH<sub>2</sub> 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](#).



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)