

# Emilin-1 monoclonal antibody (#C11A8)

## Ordering Information

[Order Online »](#)

RLT-103-M80	100µg
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	IHC (FS), WB
<b>Clone</b>	#C11A8
<b>Formulation</b>	Lyophilized
<b>Host</b>	Rat
<b>Immunogen</b>	recombinant mouse Emilin-1
<b>Reconstitution</b>	Centrifuge vial prior to opening. Reconstitute in sterile water to a concentration of 0.1-1.0 mg/ml.
<b>Species Reactivity</b>	Mouse
<b>Technical Info / Product Notes</b>	For the Original Manufacturer's data sheet please <a href="#">click here</a> .
<b>UniProt ID</b>	Q99K41



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)