

Prox1 (human), pAb

Ordering Information

[Order Online »](#)

RLT-102-PA32	200µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	When reconstituted in sterile water to a concentration of >0.5 mg/ml, the antibody is stable for at least six weeks at 2-4°C in the presence of a preservative.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application	ICC, IHC (FS), WB
Application Notes	does not work in paraffin sections
Formulation	Lyophilized from PBS, pH 7.4. Contains no preservatives.
Host	Rabbit
Immunogen	Recombinant human C-terminal part (homeo domain and prospero domain) of Prox1.
Isotype	IgG
Purity Detail	Protein A purified.
Reconstitution	Reconstitute with sterile water (>0.5mg/ml).
Species Reactivity	Human, Mouse, Rat
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



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-fr@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