

Notch1 Recombinant monoclonal antibody (E6) (Mouse IgG1 λ)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Neurogenic locus notch homolog protein 1, TAN-1, HN1
Application	ELISA, IHC
Clone	E6
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	NRR (negative regulatory region) of mouse Notch1.
Isotype	IgG1 λ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human, Mouse
Specificity	Recognizes both mouse mNRR1 and human hNRR1 but not NRR2.
Technical Info / Product Notes	This is a chimeric antibody created for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P46531 (human), Q01705 (mouse)

Last modified: May 29, 2024



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)