

Notch2 Recombinant monoclonal antibody (B6) (Rabbit IgGκ)

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

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS650-0200 | 200µg |
|-----------------|-------|

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 2, AGS2, hN2 |
| Application | ELISA, IHC |
| Clone | B6 |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Rabbit |
| Immunogen | NRR (negative regulatory region) of mouse Notch2. |
| Isotype | IgGκ |
| Purity Detail | Protein A affinity purified. |
| Species Reactivity | Human, Mouse |
| Specificity | Recognizes both mouse mNRR2 and human hNRR2, but not NRR1. |

Technical Info / Product Notes This chimeric rabbit antibody was made using the variable domain sequences of the original human scFv format, for improved compatibility with existing reagents, assays and techniques.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
Q04724 (human), O35546 (mouse)
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
5546 (mouse)
Phone: +32 3 466 0420
info-be@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

Worry-free Guarantee

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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.