

# Human Ephrin Type A receptor 2 Recombinant monoclonal antibody (1C1) (Rabbit IgGκ)

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	EphA2, Epithelial cell kinase, Tyrosine-protein kinase receptor ECK
<b>Application</b>	ELISA, Flow Cytometry, IF, WB
<b>Clone</b>	1C1
<b>Formulation</b>	Liquid. In PBS containing 0.02% Proclin 300.
<b>Host</b>	Rabbit
<b>Immunogen</b>	1C1 was isolated after screening of a non-immunized human Fab phage library. Selection was based on the ability to bind soluble human recombinant EphA2 (residues 183-510). The Fab was subsequently converted to full length human IgG1.
<b>Isotype</b>	IgGk
<b>Purity Detail</b>	Protein A affinity purified.
<b>Species Reactivity</b>	Human
<b>Specificity</b>	Recognizes the N-terminal LBD (ligand-binding domain) region of the EphA2 ectodomain (IgG1 KD ~0.8 nM; Fab KD ~140 nM) and not to other related human Eph receptors such as EphA1, EphA2, and EphA4, or any EphB receptor family members.

## Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original Human IgG1 format, for improved compatibility with existing reagents, assays and techniques.

## UniProt ID

P29317

## Worry-free Guarantee

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

**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)