

CD202B (human), (recombinant) (Fc-tag)

CD202B, also known as TIE-2, is an endothelial cell-specific receptor tyrosine kinase (RTK) that is known as a functioning molecule of vascular endothelial cells. TEK comprises a subfamily of RTK with TIE, and these two receptors play critical roles in vascular maturation, maintenance of integrity and remodeling. TEK signaling is necessary for the development of the embryonic vasculature and suggests that TEK signaling may also be required for the development of the tumor vasculature.

Ordering Information

[Order Online »](#)

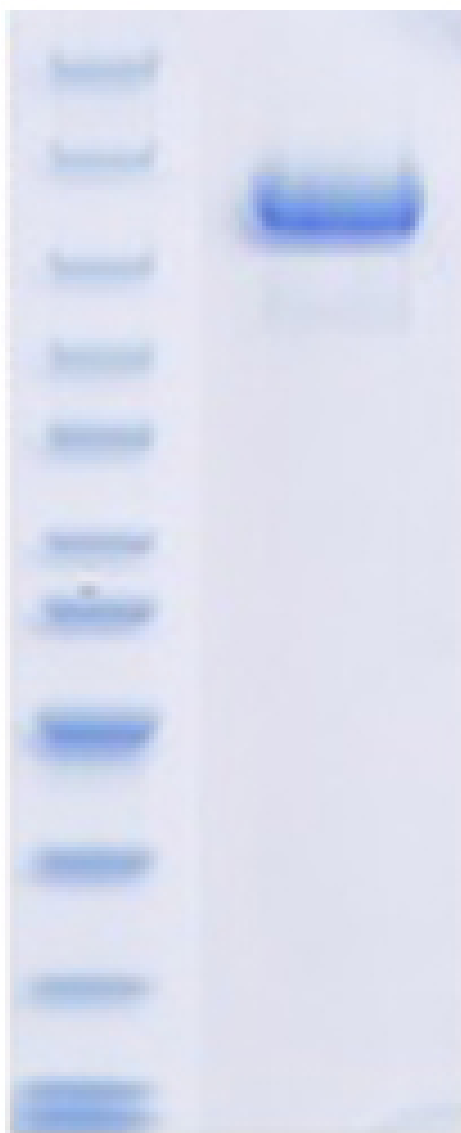
ENZ-PRT323-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa M

200
150
120
100
85
70
60
50
40
30
25
20



As a result of glycosylation, the recombinant protein migrates as an approximately 122.2 kDa and 59.5 kDa protein in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TIE-2, TEK tyrosine kinase, Angiopoietin-1 receptor

Endotoxin Content <1 EU/μg (LAL method)

Formulation Lyophilized from sterile PBS, pH 7.4.

MW ~132.5 kDa (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in HEK293 cells. A DNA sequence encoding the human CD202B (Met1-Lys745) was expressed with the fused Fc region of human IgG1 at the C-terminus.

UniProt ID Q02763



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)