

VEGFR2 (human), (recombinant) (GST- tag)

Ordering Information

[Order Online »](#)

BML-SE404-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.
Handling	After opening, prepare aliquots, freeze in liquid nitrogen and store at -80°C. Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	KDR, FLK-1, Vascular endothelial growth factor receptor 2
Application Notes	Useful tool to study enzyme regulation and kinetics, phosphorylate target substrates, screen for inhibitors.
Formulation	Liquid. In 40 mM TRIS, pH 8.0, containing 110 mM NaCl, 2.2 mM KCl, 0.02% Tween-10, 20% glycerol, and 3 mM DTT.
GenBank ID	AF035121
MW	~87 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in Sf9 insect cells with N-terminal GST tage. Human VEGFR2 (aa 805-1356). Produced in a baculovirus expression system.
Specific Activity	≥400 pmol/min/μg. Assay was done in Kinase buffer containing 0.2 mM DTT using Poly-(Glu4:Tyr)-biotin substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 40 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).
UniProt ID	P35968



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-
fr@enzolifesciences.com](mailto:info-fr@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)