VEGFR2 (human), (recombinant) (GSTtag)

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 NotesUseful 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com