

Data Sheet

Flt1 (VEGFR1)

Human, recombinant, N-terminal His tag

Catalog #: 40223

Lot#: 130121-a5 **Conc.:** 0.093 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 500 mM Imidazole, 0.04% Tween-20, 3 mM DTT, and 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Rahimi, N., *et al. J. Biol. Chem.* **275:** 16986-16992, 2000.
2. Shibuya, M. *Int. J. Biochem. Cell Biol.* **33:** 409-420, 2001.

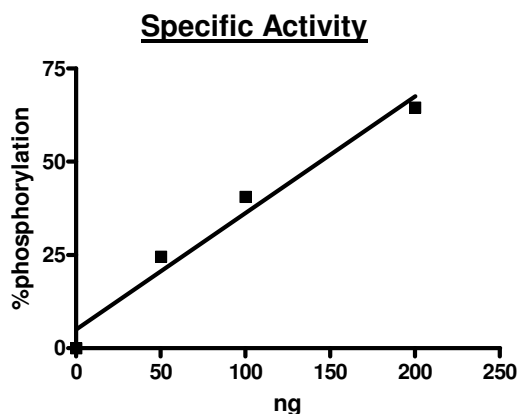
Description: Human recombinant Flt1 (FMS-like tyrosine kinase), also known as VEGFR1 (Vascular endothelial growth factor receptor 1), GenBank Accession No. AF063657, a.a. 783-end with N-terminal His tag, MW=67.5 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 7 pmol/min/μg

Assay conditions: Enzyme activity was measured using Z-LYTE assay conditions (Invitrogen). Assay buffer contains 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35, 2 μM substrate (Tyr 04, Invitrogen), and Flt1. Reaction was done at room temperature for 1 hour at 30 °C. Developer solution B was added to the reaction and incubated for 1 hour at RT before measuring the fluorescence intensity.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

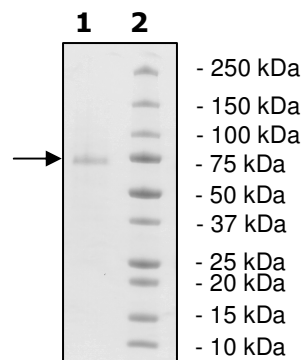
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
0.9 μg FLT1
Lane 2:
Protein Marker

MW: 67.5 kDa
Purity: ≥75%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com