

Data Sheet

FLT3

Human, recombinant, N-terminal HIS tag

Catalog #: 40225

Lot#: 1004

Conc.: 0.5 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.025% Tween-20, 125 mM Imidazole, 50% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Takahashi, S., *et al.*, *Leuk. Res.* **29 (8)**, 893-899 (2005)
2. Levis, M., *et al.*, *Blood.* **106 (2)**, 673-680 (2005).

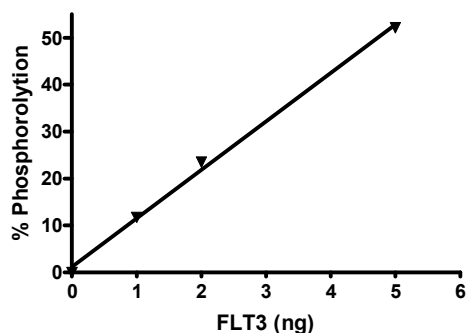
Description: Human Flt3 (GenBank Accession No. NM_004119), (a.a. 564-end) with N-terminal His tag, MW=55 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 225 pmol/min/μg.
 Assay condition: The enzyme reaction was carried out for 1h at room temperature in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EDTA, 0.01% BRIJ-35 and 200 μM of ATP. Substrate : 2 μM Tyr peptide 2 from Invitrogen.

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

Quality Assurance

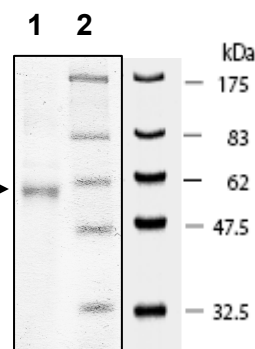
FLT3 Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
3 μg FLT3

Lane 2:
Protein Marker
BioLabs (#P7708L)



MW: 55 kDa.

Purity: ≥80%

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