

Data Sheet

c-Kit, His-tag

Human, recombinant, N-terminal His-tag

Catalog #: 40250

Lot#: 110405

Conc.: 1.3 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole, 20% Glycerol and 3 mM DTT.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Picardo M Cardinali G. J Invest Dermatol.2011 Jun;**131(6)**:1182-5.
2. Yarden, Y., *et al.* *EMBO J.* **6 (11)**,

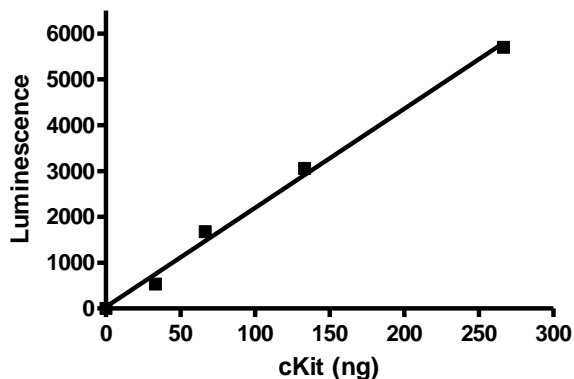
Description: Human c-Kit also known as KIT, PBT, CD117, Piebald Trait, SCFR, and Mast/Stem Cell Growth Factor Receptor, GenBank Accession No. NM_000222, a.a. 544-976(end) with N-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=53 kDa.

Specific Activity: 15 pmole/min/ μg .
 Assay Conditions: 40 mM Tris-HCl pH 7.4, 20 mM MgCl_2 , 0.1 mg/mL BSA and 1 mM DTT using 0.2 mg/ml Poly-(Glu4:Tyr)-biotin substrate and 100 μM ATP.
 Reaction was done at 30°C for 30 min.
 ATP reduction was monitored using ADP-Glo from Promega.

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

Quality Assurance

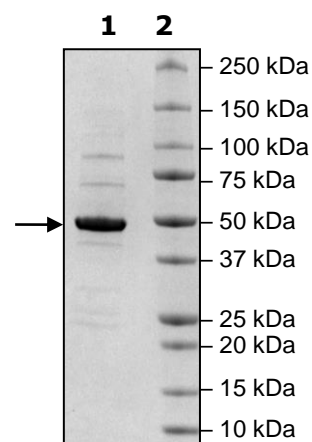
Specific Activity



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

Lane 1:
2 μg cKit
Lane 2:
Protein Marker

MW: 53 kDa
Purity: 83%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com