



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

cKit (V560G)

Human, recombinant, N-terminal His-tag

Catalog #: 40253

Lot#: 100819

Conc.: 0.05 mg/ml

Formulated in:

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

Stability:

>6 months at -80°C

References:

1. Schumaker, J.A., *et al.* (2008) *J. Clin. Pathol.* **61**:109-114
2. Sun, J. *et al.* (2009) *J. Biol. Chem.* **284**: 11039-11047.

Description: Human cKit (V560G) (GenBank Accession No. NM_000222), (a.a. 544-end) with N-terminal His tag, MW=53.4 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 100 pmol/min/μg. Enzyme reaction is conducted in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35 and 2 μM substrate (Tyr Peptide 6, Invitrogen) at room temperature for 1 hour. Purity ≥ 50%.

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

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