

## Data Sheet

### cKit (D816V), His-Tag

Human, recombinant, N-terminal His-tag

**Catalog #:** 40252

**Lot#:** 110831

**Conc.:** 0.28 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles.

#### 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.

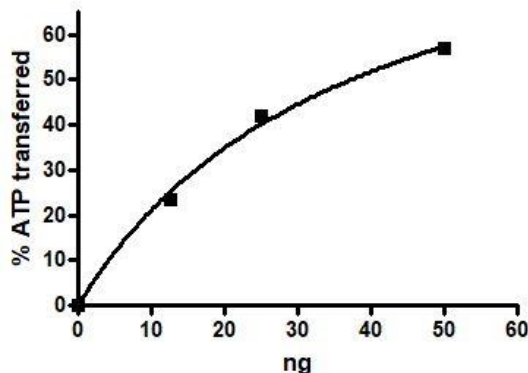
**Description:** Human cKit with D816V substitution, GenBank Accession No. NM\_000222, a.a. 544-end with N-terminal His- tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=53 kDa.

**Specific Activity:** 473 pmol/min/ $\mu\text{g}$ .  
Assay conditions: 40 mM Tris-HCl pH 7.4,  
20 mM  $\text{MgCl}_2$ , 0.1 mg/mL BSA and 1 mM  
DTT using 0.2 mg/ml Poly-(Glu4:Tyr)-biotin  
substrate and 20  $\mu\text{M}$  ATP. Reaction was  
done at  $30^{\circ}\text{C}$  for 30 min. Amount of ATP  
transferred was calculated using Kinase-Glo  
reagent from Promega

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

## Quality Assurance

### Specific Activity:



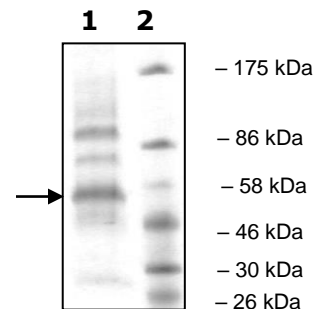
### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
4.2  $\mu\text{g}$  cKit(D816V)

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 53 kDa

**Purity:**  $\geq 32\%$



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)