

## Data Sheet

### cKit (V654A)

Human, recombinant, N-terminal His-tag

**Catalog #:** 40251

**Lot #:** 120806-G      **Conc.:** 0.27 mg/ml

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

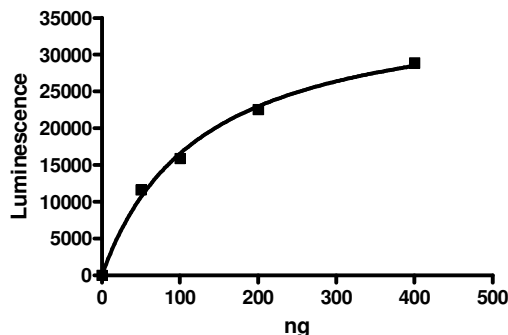
**Description:** Human Kit (V654A) (GenBank Accession No. NM\_000222), (a.a. 544-976) with N-terminal His tag, MW=53.4 kDa, expressed in an Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 89 pmole/min/μg  
 Assay buffer: 50 mM HEPES, pH 7.4, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 1 mM DTT, 3 μM Naorthovanadate, 0.1 mM ATP, 30 μg/ml Poly (Glu:Tyr)4:1 substrate, and 4 μg/ml recombinant cKit. Assay was done with 0.2 mg/ml Poly-(Glu-Tyr) peptide using ADP-Glo® (Promega) conditions at 30 °C for 35 minutes.

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

### Quality Assurance

**Specific Activity**



**10% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
 1.5 μg cKit V654A  
**Lane 2:**  
 Protein Marker  
 Biorad (#161-0373)

**MW:** 53.4 kDa  
**Purity:** ≥82%

