

## Data Sheet

### MEK1 (K97R), His-Tag

Human, Recombinant, C-terminal His-tag

**Catalog #:** 40075

**Lot#:** 190520-A

**Conc.:** 0.36 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 20% glycerol,  
3mM DTT, 140 mM Imidazole

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

**References:**

1. Smith, C.K. and Windsor, W.T.  
*Biochemistry* **46** (5), 1358-1367  
(2007).
2. Burgermeister, E. *et al. Mol. Cell. Biol.*  
**27** (3), 803-817 (2007).

**Description:** Human MEK1 (K97R) also known as Mitogen-Activated Protein Kinase Kinase 1, MAPK/ERK Kinase 1, ERK Activator Kinase 1, MAPKK 1, and PRKMK, GenBank Accession No. NM\_002755, a.a. 1-393(full length) with C-terminal His-tag, expressed in an *E. coli* expression system. MW=44 kDa.

### **Quality Assurance**

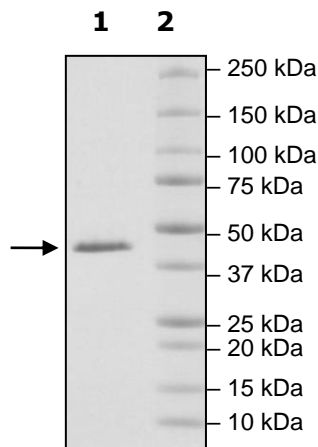
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
MEK1(K97R)

**Lane 2:**  
Protein Marker

**MW:** 44 kDa

**Purity:**  $\geq 90\%$



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 **1.858.259.5734**

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)