

Data Sheet

CSK, GST-Th-His-Tag

Human, recombinant,
N-terminal GST-Th-His-tag

Catalog #: 40410

Lot #: 1001 **Conc.:** 0.30 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 30% glycerol, 10 mM reduced glutathione, and 3 mM DTT.

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

Description: Human CSK also known as C-Terminal Src Kinase, Protein-Tyrosine Kinase CYL, C-Src Kinase and C-Src Tyrosine Kinase, GenBank Accession No. NM_004383, a.a. 1-450(full length) with N-terminal GST-Th-His-tag, expressed in Sf9 cells via a baculovirus expression system. MW = 80 kDa.

References:

1. Adachi, S. *et al.*, *Cancer Res.* **67 (13)**, 6493-6501 (2007)
2. Rosano, L. *et al.*, *Cancer Res.* **67 (13)**, 6351-6359 (2007)

Application Reference(s):

1. Figueroa, M., Oberlies, N.H., *et al.*, *The Journal of Antibiotics.* **65**, 559-564 (2012).

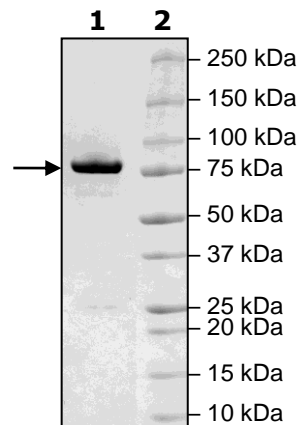
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg CSK

Lane 2:
Protein Marker

MW: 80 kDa
Purity: >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 **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com