



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

cMet (human, recombinant, N-terminal GST Tag)
Catalog #: 40255

Description: Human cMet (GenBank Accession No. X54559), (a.a. 956-1390) with N-terminal GST tag, MW= 74 kDa, expressed in an Baculovirus infected Sf9 cell expression system.

Source: Baculovirus infected Sf9 cells

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EGTA, 0.1mM EDTA, 0.1 mM PMSF, 25% glycerol.

Stability: >12 months at -70°C

Purity: > 90% by densitometry

Specific Activity: The specific activity of cMet was determined to be 46 pmol /min/ug as per activity assay protocol. One unit is defined as the amount of enzyme that will transfer 1 pmol phosphate to Tyr substrate per minute at pH 7.4 and 30°C. Assay buffer: 50 mM HEPES, pH 7.4, 3 mM MgCl₂, 3 mM MnCl₂, 1 mM DTT, 3 uM Na-orthovanadate, 0.1 mM ATP, 30 ug/ml Poly (Glu:Tyr)4:1 substrate, and 4 ug/ml recombinant cMet

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

References:

1. Giordano, S. et al., Oncogene, 1989, 4: 1383-1388.
2. Lyer, A. et al., Cell Growth Differ. 1990, 1: 87-95.

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** or **1.858.829.3082**, Fax **1.858.259.5734**

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

Please visit our website at: www.bpsbioscience.com