



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

EphB4, active

(Recombinant human protein expressed in Sf9 cells)

Catalog # **E-K062**
Size: 10 µg
Specific activity: 125 U/mg

Product Description

EphB4 (EPH receptor B4) is a receptor tyrosine kinase. Recombinant human EphB4 (amino acids 563-987) with N-terminal 6xHis tag was expressed in a baculovirus infected Sf9 insect cell expression system. The GenBank Accession number is NM_004444. MW = 50 kDa.

Specific Activity

125 U/mg. One unit of EphB4 activity is defined as 1 nmol phosphate incorporated into tyrosine substrate per minute at pH 7.4 and 30°C.

Assay Condition

50 mM HEPES, pH 7.4, 3 mM MgCl₂, 3 mM MnCl₂, 1 mM DTT, 3 µM Na-orthovanadate, 0.1 mM ATP, 30 µg/ml Poly (Glu:Tyr) 4:1 substrate, and 0.4 µg/ml recombinant EphB4.

Purity

> 90% by SDS-PAGE.

Formulation

Recombinant protein is in storage buffer (25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 3 mM DTT, and 50% glycerol).

Storage and Stability

Store product frozen at or below -70 °C. Enzyme will be stable for at least 6 months at -70 °C as undiluted stock. Freeze in working aliquots to avoid repeated thawing and freezing.

Application

EphB4 is useful for the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

References

1. Bruhl T. *et al.* (2004) *Circ. Res.* **94** (6), 743-751
2. Fuller T. *et al.* (2003) *J. Cell. Sci.* **116** (Pt 12), 2461-2470