



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.

EphA2, active

(Recombinant human protein expressed in Sf9 cells)

Catalog # **E-K060**
Size: 10 µg
Specific activity: 170 U/mg

Product Description

EphA2 (EPH receptor A2) is a receptor tyrosine kinase. Recombinant human EphA2 (amino acids 596-end) with N-terminal 6xHis tag was expressed in a baculovirus infected Sf9 insect cell expression system. The GenBank Accession number is NM_004431. MW = 47 kDa.

Specific Activity

170 U/mg. One unit of EphA2 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 0.1 µg/ml recombinant EphA2.

Purity

> 80% 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

EphA2 is useful for the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

References

1. Macrae M. *et al.* (2005) *Cancer Cell* **8** (2), 111-118
2. Xu H. *et al.* (2005) *Am. J. Physiol. Renal Physiol.* **288** (4), F855-F866
3. Walker-Daniels J. *et al.* (2003) *Am J Pathol*, **162**, (4), 1037-1042

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

TDS E-K060 Rev: 1 (03/15/06)