

Data Sheet

GSK3 β , GST-tag

Human, recombinant, N-terminal GST-tag

Catalog #: 40007

Lot #: 150501-G2

Conc.: 0.18 mg/ml

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

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

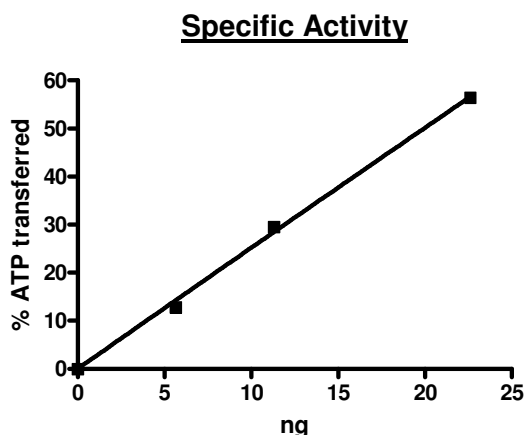
1. Sutherland C. *et al. Biochem J.* 1993 Nov 15; **296** (Pt 1):15-9.
2. Sperber BR *et al. Neurosci Lett.* 1995 Sep 8; **197**(2):149-53.

Description: Human GSK3 β (GenBank Accession No. NM_002093), full length, with N-terminal GST tag, MW=73 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: ≥ 500 pmol/min/ug. Enzyme reactions include 50 μL volumes of 1x Kinase buffer (40 mM Tris pH 7.4, 20 mM MgCl_2 , 0.1 mg/mL BSA) along with 20 μM ATP, 0.2 mg/mL GSK-tide substrate, 3mM DTT, and GST-GSK3 β . After 45 min incubation at 30°C , 50 μL of the Kinase-Glo reagent solution was added and luminescence from remaining ATP was measured.

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

Quality Assurance

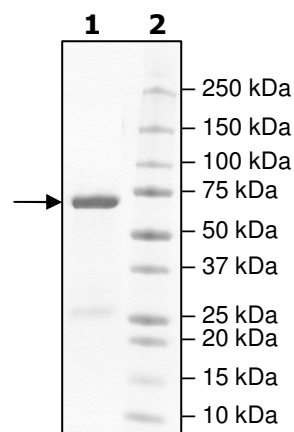


4-20% SDS-PAGE Coomassie staining

Lane 1:
2.5 μg GSK3 beta

Lane 2:
Protein Marker

MW: 74.5 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com