

Data Sheet

GSK3 α , GST-tag

Human, recombinant, N-terminal GST tag

Catalog #: 40006

Lot #: 130613-H1 **Conc.:** 1.0 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 4 mM GSH, 3 mM DTT, 20% glycerol

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

References:

1. Meijer, L. *et al.*, *Trends Pharmacol Sci.* **25 (9)**, 471-480 (2004).
2. Patel, S. *et al.*, *Biochem Soc. Trans.* **32(pt5)**: 803-808 (2004).

Description: Human GSK3 α , also known as GSK3 α , glycogen synthase kinase 3 alpha, GSK3alpha, GenBank Accession No. BC027984, a.a. 2-483(end), with an N-terminal GST-tag, expressed in a Baculovirus infected Sf9 cell expression system, MW=78 kDa.

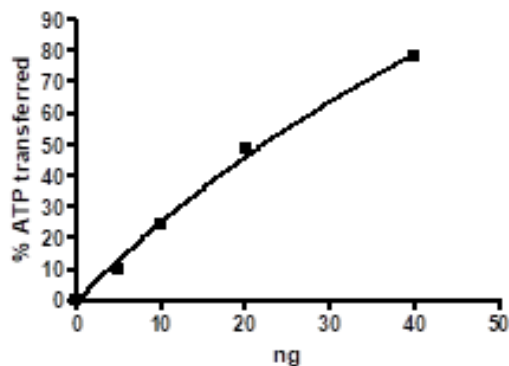
Specific Activity: ≥ 500 pmol/min/ug.

Assay Conditions: Enzyme reactions include 50 μL volumes of 1x Kinase buffer (40 mM Tris pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA) along with 20 μM ATP, 0.2 mg/mL GSKtide substrate, 3mM DTT, and GST-GSK3 α . After 45 min incubation at 30 $^{\circ}\text{C}$, 50 μL of the Kinase-Glo reagent (Promega) 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

Specific Activity

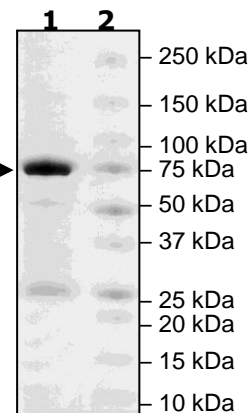


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg GSK3 α

Lane 2:
Protein Marker

MW: 78 kDa
Purity: $\geq 81\%$



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