

Data Sheet

GSK α , GST-Tag, His-tag

Human, Recombinant, N-terminal GST-tag,
 C-terminal His-tag

Catalog #: 100814

Lot #: 200729-G2 **Conc.:** 0.56 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 3 mM DTT, and 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).

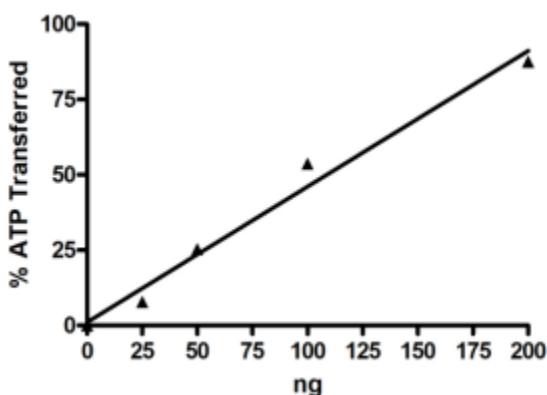
Description: Human glycogen synthase kinase 3 alpha (GSK3 α), also known as GSK3a and GSK3alpha, GenBank Accession No. NM_019884, a.a. 2-483(end), with N-terminal GST-tag and C-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system, MW=78 kDa.

Specific Activity: 109 pmol/min/ μg
 Assay Conditions: Assay was performed according to the GSK3 α Assay Kit (BPS Bioscience, #79698) with GSK3a titrated from 200 ng/rxn– 0 ng/rxn.

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

Quality Assurance

Specific Activity

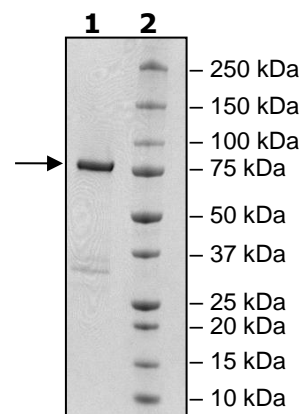


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg GSK α
Lane 2:
Protein Marker

MW: 78 kDa

Purity: 85%



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** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com