

GSK3-beta

glycogen synthase kinase 3 beta

Recombinant Human Active Protein Kinase

HGNC Symbol: GSK3B

Synonyms: n/a

Product No.: 0310-0000-1

Lot: 003

Description: Human GSK3-beta, full-length, amino acids M₁-T₄₂₀ (as in [NCBI/Protein](#) entry NP_001139628.1), N-terminal GST-HIS₆ fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: GSK3-beta Lot 003, was confirmed as GSK3-beta by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 76,312 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.276 µg/µl

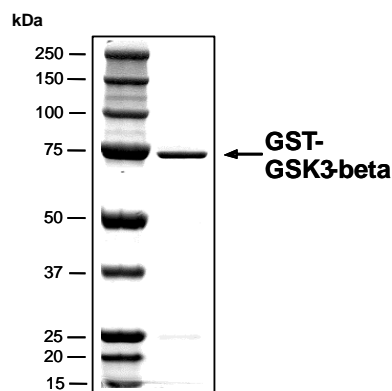
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 6.3 pmol/µg × min

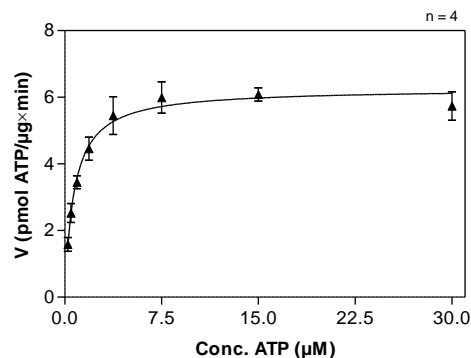
ATP-K_M: 0.7 µM

**GSK3-beta Lot 003:
Coomassie stain**



2.0 µg GST-GSK3-beta

**GSK3-beta Lot 003:
Determination of V_{max} and K_M value for ATP**



Determination of K_M value & Specific activity:

- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg/ml PEG_{20,000}
 - ATP (variable)
 - Substrate: RBER-IRStide 100 µg/ml
 - Kinase: 2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

GSK3-beta

Product No.: 0310-0000-1

GST-GSK3-beta Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG HHHHHG	RRRASVAAGI	240
241	LVPRGS PGLD	GIYARE MSGR	PRTTSFAESC	KPVQPSAFG	SMKVS RDKD G	SKVTTV VATP	300
301	GQGPDR PQEV	SYTDTKVIGN	GSFGVVYQAK	LCDSGELVAI	KKVLQDKRFK	NRELQIMRKL	360
361	DHCNIV RLRY	FFYSSGEK KD	EVYLN LVL LDY	VPETVYR VAR	HYSRAKQ TL P	VIYVKLY MYQ	420
421	LFRSLAY IHS	FGICHR DIK P	QNL LLDP DTA	VLKLCDF GSA	KQLVRGE PNV	SYICSR YYRA	480
481	PELIFG ATDY	TSSIDV WSAG	CVLAELL L GQ	PIFPGD SGVD	QLVEI IKVL G	TPTREQ I REM	540
541	NPNYTE FKFP	QIKAHP WTKV	FRPRT PPEAI	ALCSRL LEYT	PTARLT PLEA	CAHSFF DEL R	600
601	DPNVK LPNGR	DTPALF NFTT	QELSS NPPLA	TILIP PHARI	QAAAS TPTNA	TAASD ANT GD	660
661	RGQT NAASA	SAS NST					720

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **blue**: GSK3-beta

GSK3-beta wt ¹ Amino Acid Sequence							
1	MSGR PRTTSF	AESCK PVQQP	SAFG SMK VSR	DKDG SKV TV	VATP GQGPDR	PQEV SYTD TK	60
61	VIGN SFGVV	YQAK LCDSGE	LVAI KKVL QD	KRFK NREL QI	MRKL DHCNIV	RLRY FF YSSG	120
121	EKKD EVY LNL	VLDY VPETVY	RVAR HYSRAK	QTL PVIY VKL	YMYQ LFRSLA	YIHS FGI CHR	180
181	DIK QNL LLD	PDTAV LK LCD	FGSA Q LV RG	EPNV SYI CSR	YYRA PEL IFG	ATDY TSS IDV	240
241	WSAG CVL AEL	LLGQ PI FPGD	SGVD Q LVEI	KVLG T P TREQ	I REMNP NYTE	FKFP Q IKAHP	300
301	WTKV FR PRTP	PEAI ALCSRL	LEYT P TARLT	PLEA CAHS FF	DEL RDPNV KL	PNGR D T PALF	360
361	NFTT Q ELSS N	PPLAT IL IP	HARI Q AAAS T	PTNA T AASDA	NTGD R GQ TNN	AASAS AS NS T	420

blue: GSK3-beta sequence expressed in recombinant protein

¹[NCBI/Protein](https://www.ncbi.nlm.nih.gov/Protein) accession number NP_001139628.1

This product was manufactured at ProQinase in Freiburg, Germany, and is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.