

ProQinase™ BMPR1B

bone morphogenetic protein receptor type 1B

Recombinant Human Active Protein Kinase

HGNC Symbol: BMPR1B

Synonyms: ALK6, CDw293

Product No.: 1667-0000-1

Lot: 001

Description: Human BMPR1B, C-terminal fragment, amino acids Y₁₄₉-L₅₀₂ (as in [NCBI/Protein](#) entry NP_001194.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

Product identity: BMPR1B Lot 001, was confirmed as BMPR1B by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 63,387 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM HEPES pH 7.5, 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: 0.674 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

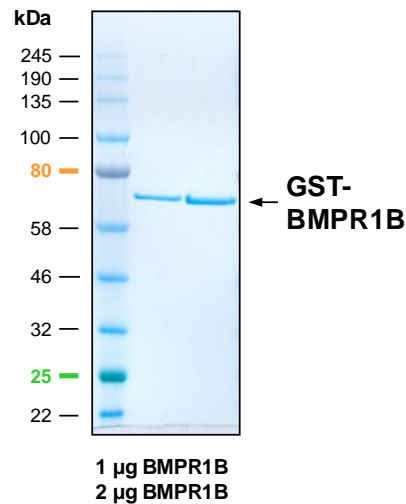
Biochemical Parameters:

Specific kinase activity (P_i transfer): 13 pmol/µg × min
ATP-K_M: 0.2 µM

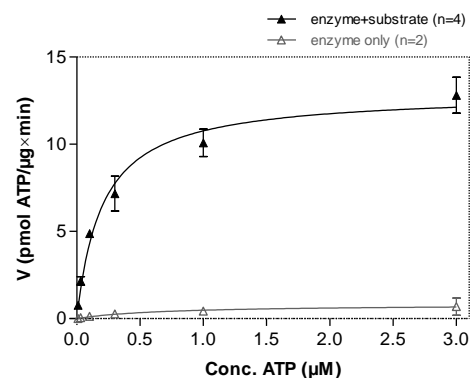
Additional assay technology:

BMPR1B Lot 001 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

BMPR1B Lot 001: Coomassie stain



BMPR1B Lot 001: 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: Casein 40 µg/ml
 - Kinase: 0.2 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

ProQinase™ BMPR1B

Product No.: 1667-0000-1

GST-BMPR1B 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	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVL FQG	240
241	PKRQETRPR	YSIGLEQDET	YIPPGESLRD	LIEQSQSSGS	GSGPLLLVQR	TIAKQIQMVK	300
301	QIGKGRYGEV	WMGKWRGEKV	AVKVFVFTTEE	ASWFRETEIY	QTVLMRHENI	LGFIADIKG	360
361	TGSWTQLYLI	TDYHENGSLY	DYLNKSTTLDA	KSMKLKAYSS	VSGLCHLHTE	IFSTQKPAI	420
421	AHRDLKSKNI	LVKKNGTCCI	ADLGLAVKFI	SDTNEVDIPP	NTRVGTKRYM	PPEVLDES LN	480
481	RNHFSYI MA	DMYSFGLILW	EVARRCVSGG	IVEEYQLPYH	DLVPSDPSYE	DMREIVCIKK	540
541	LRPSFPNRWS	SDECLRQMGK	LMTECWAHNP	ASRLTALRVK	KTLAKMSESQ	DIKL	600

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: BMPR1B fragment

BMPR1B wt ¹ Amino Acid Sequence							
1	MLRSAGKLN	VGTKKEDGES	TAPTPRPKVL	RCKCHHHCPE	DSVNNICSTD	GYCFTMIEED	60
61	DSGLPVVTS G	CLGLEGSDFQ	CRDTPIPHQR	RSIECCTERN	ECNKDLHPTL	PPLKNRDFVD	120
121	GPIHHRALLI	SVTVC SLLLV	LIILFCYFRY	KRQETRPRYS	IGLEQDETYI	PPGESLRDLI	180
181	EQSQSSGSGS	GLPLL VQRTI	AKQIQMVKQI	GKGRYGEVWM	GKWRGEKVAV	KVFFVTEEAS	240
241	WFRETEIYQT	VLMRHENILG	FIAADIKGTG	SWTQLYLITD	YHENGSLYDY	LKSTTLDAKS	300
301	MLKLAYSSVS	GLCHLHTEIF	STQKPAIAH	RDLKSKNILV	KKNGTCC IAD	LGLAVKFISD	360
361	TNEVDIPPNT	RVGTRKYMPP	EVLDES LN RN	HFQSYIMADM	YSFGLILWEV	ARRCVSGGIV	420
421	EEYQLPYHDL	VPSDPSYEDM	REIVCIKKLR	PSFPNRWSSD	ECLRQMGKLM	TECWAHNPAS	480
481	RLTALRVKKT	LAKMSESQDI	KL				540

blue: BMPR1B sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_001194.1