

Data Sheet

CDK5/p35

Human, recombinant, N-terminal GST tag

Catalog #: 40095

Lot #: 121101 **Conc.:** 0.63 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Tang, D., *et al.*, *Prog Cell Cycle Res.* 1996;**2** :205-16.
2. Dhavan, R., *et al.*, *Nat Rev Mol Cell Biol.* 2001 Oct;**2**(10):749-59.

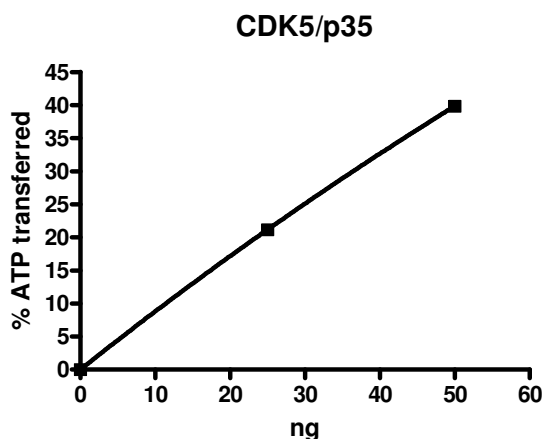
Description: Complex of N-terminal GST-tagged recombinant full length human CDK5 (GenBank Accession No. NM_004935), and N-terminal GST-tagged recombinant full length human p35 (GenBank Accession No. NM_003885). Co-expressed in a Baculovirus-infected Sf9 cell expression system. CDK5 MW=60 kDa, p35 MW=61 kDa.

Specific Activity: 180 pmol/min/μg
 Assay conditions: Assay was done in Kinase buffer containing 0.2 mM DTT using Histone 1 substrate (0.2 mg/ml) and 20 mM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo[®] reagent (Promega).

Application: Useful for three-dimensional structure studies.

Quality Assurance

Specific Activity



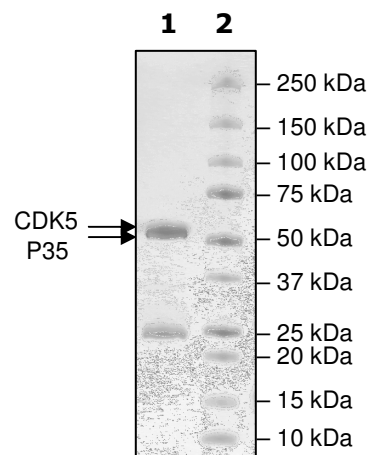
4-20% SDS-PAGE Coomassie staining

Lane 1:
1.8 μg CDK5/p35

Lane 2:
Protein Marker

MW:
CDK5: 60 kDa
p35: 61 kDa

Purity: ≥60%



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

CDK5

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDPLVPRGSPGFMQKY
EKLEKIGEGTYGTVFKAKNRETHEIVALKRVRLLDDDEGVPSSALREICLLKELKHKNI
VRLHDVLHSDKCLTLVFEFCDQDLKKYFDFSCNGDLDPKIVKSLFQLLKGLGFCHSRN
VLHRDLKPNLLINRNGELKLADFGARAFGIPVRCYSAEVVTLWYRPPDVLFGAKLY
STSIDMWSAGCIFAELANAGRPLFPGNDVDDQLKRIFRLLGTPTEEQWPSMTKLPDY
KPYPMPATTSLVNVVVKLNATGRDLLQNLKCNPVQRISAEALQHPYFSDFCPP

P35

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDPLVPRGSPGFMGTV
LSLSPSYRKATLFEDGAATVGHYTAQNSKNAKDKNLKRHSISVLPWKRIVAVSAKK
KNSKKVQPNSSYQNNITHLNNENLKKSLSCANLSTFAQPPPAQPPAPPASQLSGSQT
GGSSSVKKAPHPAVTSAGTPKRVIVQASTSELLRCLGEFLCRRCYRLKHLSPDTPVL
WLRSDRSLLLQGWQDQGFITANVVFLYMLCRDVISSEVGSDELQAVLLTCLYLS
YSYMGNEISYPLKPLVESCKEAFWDRCLSVINLMSSKMLQINADPHYFTQVFSDLKN
ESGQEDKKRLLLGLDR

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