

Product Information

LRRK2[G2019S]

Product Number : **09-146**

Product description

Human LRRK2, catalytic domain [960-2527(end) amino acids and G2019S of accession number NP_940980.3] was expressed as N-terminal GST-fusion protein (206 kDa) using baculovirus expression system. GST-LRRK2[G2019S] was purified by using glutathione sepharose chromatography and gel filtration chromatography.

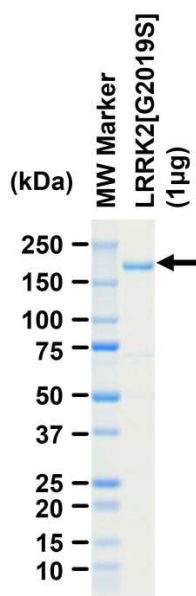
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
Avoid repeating freeze-thaws.

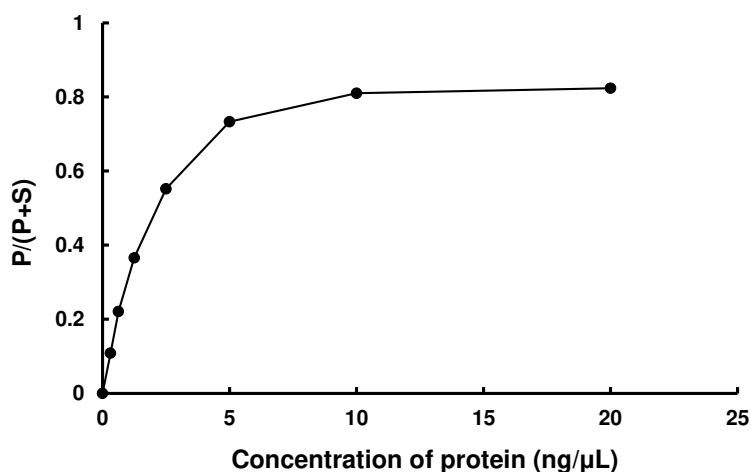
SDS-PAGE



Purity: 94 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : LRRKtide

ATP : 500 μM