

Product Information

LRRK2

Product Number : **09-110**

Product description

Human LRRK2, catalytic domain [960-2527(end) amino acids of accession number AAV63975.1] was expressed as N-terminal GST-fusion protein (206 kDa) using baculovirus expression system. GST-LRRK2 was purified by using glutathione sepharose chromatography.

*LRRK2(09-110) is not available outside Japan.

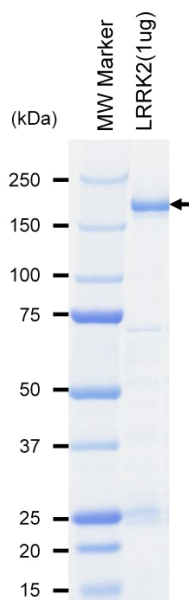
Storage buffer:

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.

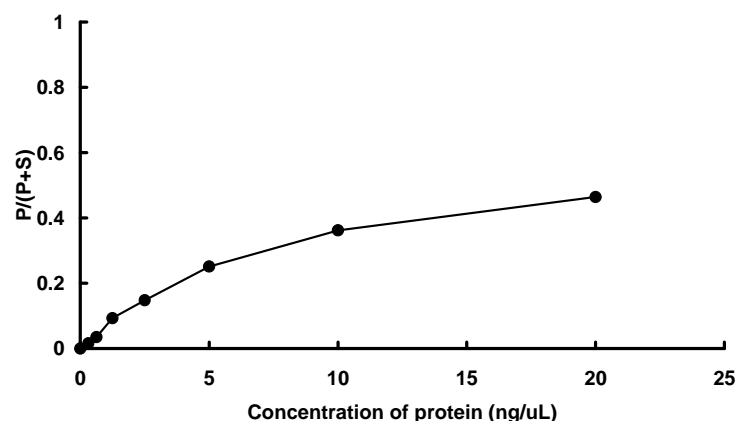
SDS-PAGE



Purity: 56 %

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 μ M