

Data Sheet

Cas9, His-tag (*S. pyogenes*)

Human recombinant, N-terminal His-tag

Catalog #: 100206

Lot #: 181025-G **Conc:** 1.46 mg/ml

Formulated in: 20 mM HEPES pH 7.5, 500 mM KCl, 20% Glycerol, 3 mM DTT.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Liang, X. *et al.*, *J Biotechnol.* 2017; **241**:136-146
2. Kim, S., *et al.*, *Genome Res.* 2014; **24(6)**: 1012-1019

Description:

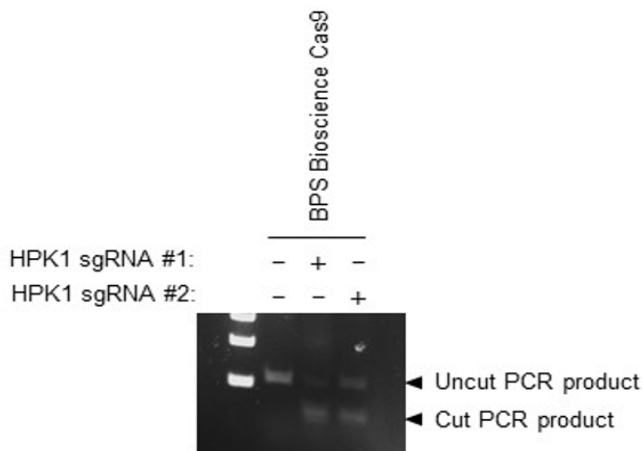
Streptococcus pyogenes CRISPR associated protein 9 (Cas9), also known as SpCas9, GenBank Accession No. NC_002737.2, a.a. 2-1368(end) with N-terminal His-tag, expressed in an *E. coli* cell expression system. MW = 160 kDa.

Assay Conditions:

The Cas9 nuclease was incubated with sgRNA at 37°C for 10 minutes to form a functional RNP complex. A purified PCR product was then added at 37°C for 60 minutes, before Cas9 activity was visualized by agarose gel.

Applications: Cas9 can be used to form CRISPR RNP complexes for in vitro cleavage assays to identify sgRNAs with the highest cutting efficiencies.

Quality Assurance

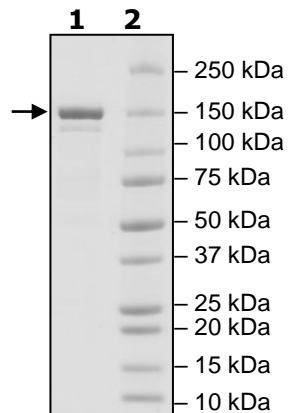


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 CAS9
Lane 2:
 Protein Marker

Purity: $\geq 90\%$

MW: 160 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com