

Cas9 (D10A, H840A), NLS, His-Tag (*S. pyogenes*) Recombinant

Catalog: 101633
Lot: 230110

Product Information

Description:	Recombinant <i>S. pyogenes</i> SpCas9(D10A, H840A)-NLS (conserved hypothetical protein), full length. This construct contains a C-terminal His-Tag. SpCas9(D10A, H840A)-NLS is a catalytic mutant of Cas9 that contains a nuclear localization sequence for optimal activity.
Species:	<i>Streptococcus pyogenes</i>
Construct:	Cas9 (D10A, H840A) (Full Length-NLS-His) (<i>S. pyogenes</i>)
Mutation:	D10A, H840A
Concentration:	0.20 mg/ml
Expression System:	<i>E. coli</i>
Purity:	80%
Format:	Aqueous buffer solution.
Formulated In:	50 mM sodium phosphate, pH 7.5, 300 mM NaCl, 1 mM DTT, 200 mM imidazole and 10% glycerol
MW:	160 kDa
Genbank Accession:	AAK33936
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

