

## Product Information

<b>Description:</b>	Recombinant <i>S. pyogenes</i> SpCas9(H840A)-NLS (conserved hypothetical protein), full length. This construct contains a C-terminal His-Tag. SpCas9(H840A)-NLS is a catalytic mutant of Cas9 that contains a nuclear localization sequence for optimal activity.
<b>Species:</b>	<i>Streptococcus pyogenes</i>
<b>Construct:</b>	Cas9 (H840A) (Full Length-NLS-His) ( <i>S. pyogenes</i> )
<b>Mutation:</b>	H840A
<b>Concentration:</b>	0.20 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	70%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM sodium phosphate, pH 7.5, 300 mM NaCl, 1 mM DTT, 200 mM imidazole and 10% glycerol
<b>MW:</b>	160 kDa
<b>Genbank Accession:</b>	AAK33936
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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

