

## Product Information

<b>Description:</b>	Recombinant <i>Salmonella typhimurium</i> sialidase, also known as NANH (N-acylneuraminase glycohydrolase), full length encompassing amino acids 2-382. The construct contains an N-terminal His-tag (6xHis). This protein was affinity purified.
<b>Background:</b>	Sialidases are glycosyl hydrolases that cleave the $\alpha$ -ketosidic bond from sialic acid residues. These proteins are crucial for microbes like Salmonella to successfully infect their hosts by degrading the cell's glycans and contributing to adherence and infection. <i>Salmonella typhimurium</i> infection results in salmonellosis, characterized by gastrointestinal symptoms, such as abdominal cramps and diarrhea, which last from 4 to 7 days. It can, however, result in severe cases requiring hospitalization. Inhibitors targeting sialidases may prove useful in the treatment of bacterial infections.
<b>Species:</b>	<i>Salmonella Typhimurium</i>
<b>Construct:</b>	Sialidase (His-2-382(end))
<b>Concentration:</b>	0.90 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	$\geq 90\%$
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	43 kDa
<b>Genbank Accession:</b>	AE006468.2
<b>Stability:</b>	At least 6 months at $-80^{\circ}\text{C}$ .
<b>Storage:</b>	$-80^{\circ}\text{C}$
<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

