

## Product Information

<b>Description:</b>	Recombinant NEU2 (Neuroaminidase 2 or sialidase-2), full length, encompassing amino acids 2-380(end). This construct contains an N-terminal His-Tag (6xHis). This recombinant protein was affinity purified.
<b>Background:</b>	NEU2, also known as neuroaminidase 2 or sialidase-2, is a glycohydrolytic enzyme involved in the removal of sialic acid from either proteins or lipids. This enzyme is present at low levels in most tissues but plays a crucial role in carcinogenesis. Cytosolic NEU2 regulates the PI3K-Akt/mTOR pathway and is thus involved in apoptosis, drug resistance, and cancer cell migration.
<b>Species:</b>	Human
<b>Construct:</b>	NEU2 (His-2-380(end))
<b>Concentration:</b>	0.67 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥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>	NM_005383.2
<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

