

## Product Information

<b>Description:</b>	Recombinant human NPM1 (nucleophosmin 1), full length. This construct contains an N-terminal His-tag and was affinity purified.
<b>Species:</b>	Human
<b>Construct:</b>	NPM1 (His-Full Length)
<b>Concentration:</b>	0.10 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	80%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol
<b>MW:</b>	40 kDa
<b>Genbank Accession:</b>	NM_002520
<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

