

## Product Information

<b>Description:</b>	Recombinant human NPM1 (nucleophosmin), encompassing amino acids 2-294(end). This construct contains an N-terminal His-tag (6xHis) followed by a TEV site. This protein was affinity purified.
<b>Background:</b>	NPM1 (nucleophosmin), also known as nutramin or nucleolar phosphoprotein B23, is a protein that binds to nucleolar ribonucleoprotein structures and is involved in ribosome biogenesis, centrosome duplication, RNA helix destabilization, DNA repair, and tumor suppression. The encoded phosphoprotein shuttles between the nucleolus, nucleus, and cytoplasm, chaperoning ribosomal proteins and core histones from the nucleus to the cytoplasm. This protein is also known to sequester the tumor suppressor ARF in the nucleolus, protecting it from degradation until it is needed. Abnormalities in the NMP1 gene have been found in non-Hodgkin lymphoma, acute promyelocytic leukemia and acute myeloid leukemia (AML). In AML this protein was found to be mutated, lacking a folded C-terminal domain. The development of molecules able to either fold the protein to its native state or remove it from the cytosol may be beneficial in cancer therapy.
<b>Species:</b>	Human
<b>Construct:</b>	NPM1 (His-TEV-2-294(end))
<b>Concentration:</b>	0.62 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, 500 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% glycerol, and 238 mM Imidazole
<b>MW:</b>	34 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.
<b>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

