

## Product Information

<b>Construct:</b>	CDK13 (GST-516-1512(end)) / Cyclin K (GST-2-580(end))
<b>Concentration:</b>	0.59 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
<b>Expression System:</b>	Sf9; Coexpressed with CAK1, <i>S. Cerevisiae</i> (FLAG-2-368(end))
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	CDK13: NM_003718; Cyclin K: NM_001099402
<b>MW:</b>	CDK13: 138 kDa; Cyclin K: 90 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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

