

Product Information

Construct:	Transglutaminase (29-262(end)-His) (<i>K. albida</i>)
Concentration:	0.28 mg/ml
Species:	<i>Kutzneria albida</i>
Formulated In:	40 mM Tris pH8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Uniprot:	W5WHY8
MW:	28 kDa
Purity:	>90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

