

Product Information

Description:	Recombinant <i>K. albida</i> transglutaminase, encompassing amino acids 29-262end. This construct contains a C-terminal His-tag (6xHis). The recombinant protein was affinity purified.
Species:	<i>Kutzneria Albida</i>
Construct:	Transglutaminase (29-262(end)-His) (<i>K. albida</i>)
Concentration:	0.28 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	28 kDa
Uniprot:	W5WHY8
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

