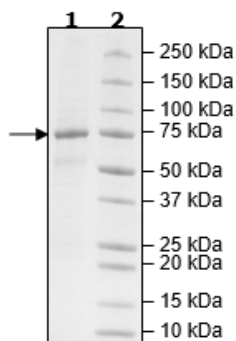


Product Information

Construct: PADI2 (1-665(end)-FLAG)
Concentration: 0.46 mg/ml
Species: Human
Formulated In: 40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 80 ng/μl FLAG peptide and 20% Glycerol.
Expression System: Sf9
Format: Aqueous buffer solution
Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage: -80°C
Genbank Accession: NM_007365
MW: 77 kDa
Purity: 84%
Specific Activity: ≥19 pmol/min/μg
Assay Conditions: Enzyme reaction was done in a buffer (50 mM HEPES, pH 7.4, 50 mM NaCl, 10 mM CaCl₂, and 2 mM DTT) containing PADI and 10 mM BAEE as a substrate for 30 min at 37°C. Amount of ammonia produced was measured and used for specific activity calculation.
 After 30 min incubation at 37°C, the reaction was quenched using 15 μL of 500 mM EDTA (final concentration is 115 mM) followed by measuring the amount of NH₃ produced by o-phthalaldehyde (OPT) and dithiothreitol (DTT). Fluorescence was measured at λ_{ex}413 nm and λ_{em}476 nm.
Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



PADI2 Activity

