

## Data Sheet

### PADI3

Human, recombinant, C-terminal FLAG-tag

**Catalog #:** 50153

**Lot#:** 110607      **Conc.:** 0.25 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/ul Flag peptide, 20% Glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

#### References:

1. Kanno T, *et al.* (2000). *J. Invest. Dermatol.* **115** (5): 813–23.
2. Nachat R, *et al.* (2005). *J. Invest. Dermatol.* **124** (2): 384–93
3. Dong S, *et al.* (2006). *Biochem. J.* **397** (3): 449–59.

#### Description:

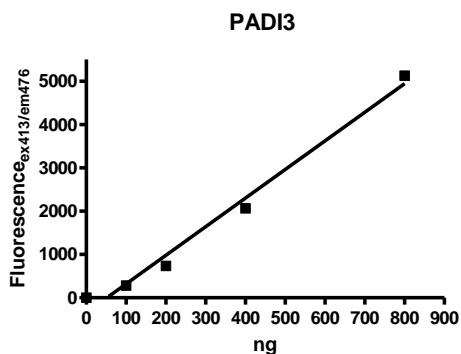
Human PADI3, also known as Peptidyl arginine deiminase, type III, (GenBank Accession No. NM\_016233), 1-664(end), with C-terminal FLAG-tag, MW= 75 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

#### Specific Activity: 22 pmole/min/μg

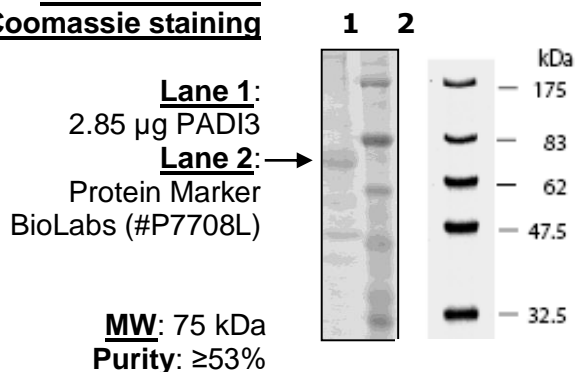
Assay Conditions: Enzyme reaction took place in a buffer (50 mM Tris, pH 7.4, 50 mM NaCl, 10 mM CaCl<sub>2</sub>, and 2 mM DTT) containing PADI and 10 mM BAEE as a substrate for 25 min at 37°C.

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 NH<sub>3</sub> produced by o-phthalaldehyde (OPT) and dithiothreitol (DTT). Fluorescence was measured at λ<sub>ex</sub>413 nm and λ<sub>em</sub>476 nm.

### Quality Assurance



#### 10% SDS-PAGE Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
 Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)