

## Data Sheet

### PADI1

Human, recombinant, C-terminal FLAG-tag

**Catalog #:** 50151

**Lot#:** 100914      **Conc.:** 0.29 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 90 µg/ml  
 FLAG peptide, 3 mM DTT and 10% glycerol.

**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. Guerrin M, *et al.* (2003). *Biochem. J.* **370** (Pt 1): 167–74.

**Description:** Human PADI1 (GenBank Accession No. NM\_013358), 1-663(end), with C-terminal FLAG-tag, MW= 75 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

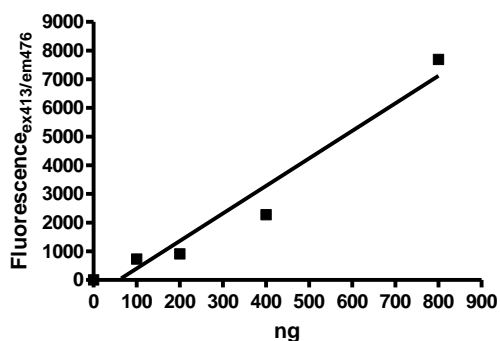
**Specific Activity:** 35 pmol/min/µg

Enzyme reaction was done in a buffer (50 mM Tris, pH 7.4, 50 mM NaCl, 10 mM CaCl<sub>2</sub>, and 2 mM DTT) containing PADI1 and 10 mM BAEE as a substrate for 25 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 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

#### Specific Activity

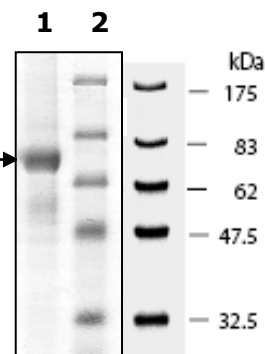


#### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
5 µg PADI1

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 75 kDa.  
**Purity:** >70%



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)