

## 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.

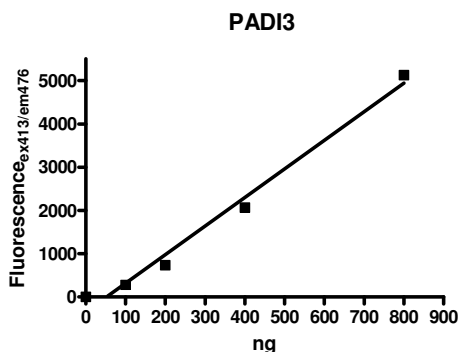
**Specific Activity:** 22 pmole/min/μg

**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.

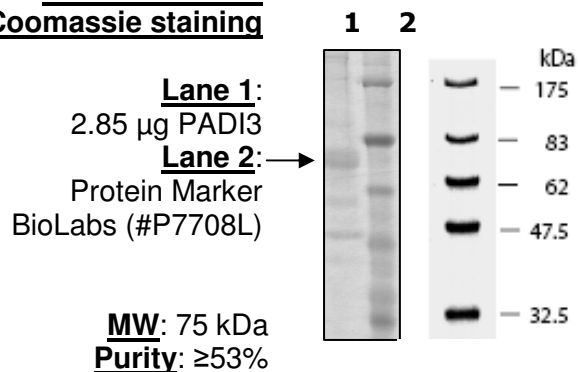
#### Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### 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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)