

Data Sheet

PADI2

Human, recombinant, C-terminal FLAG-tag

Catalog #: 50152

Lot#: 111018

Conc.: 0.06 mg/ml

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

Stability: >6 months at -80°C

References:

1. Dong S, et al. (2005.) *J. Invest. Dermatol.* 124(5):1026-33.
2. Wantanabe Y, et al. (2009). *J Hum Genet.* 54 (7): 430-2.

Description:

Human PADI2, also known as Peptidyl arginine deiminase, type II, (GenBank Accession No. NM_007365), 1-665(end), with C-terminal FLAG-tag, MW= 77 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

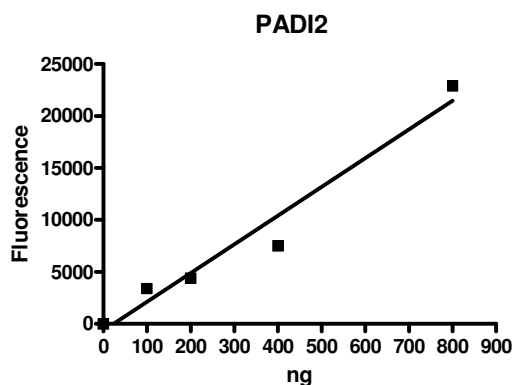
Specific Activity: 19 pmole/min/μg.

Assay Conditions: Enzyme reaction took place in a buffer (50 mM Tris, 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.

Application:

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

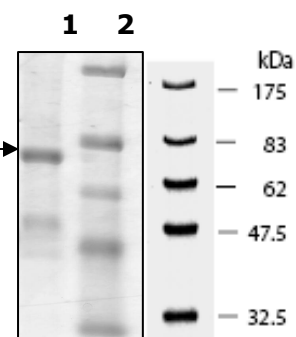
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 2.6 μg PADI2
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 77 kDa
Purity: ≥61%



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

Please visit our website at: www.bpsbioscience.com