

PAD6 Protein

Full length recombinant human protein expressed in Sf9 cells

Catalog # P312-30FG

Lot # T1065-3

Product Description

Recombinant full-length human PAD6 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_207421](#).

Gene Aliases

PADI6

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

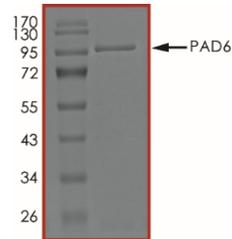
Scientific Background

PAD6 or PADI6 is a member of the peptidylarginine deiminases which convert arginine residues to citrulline residues in the presence of calcium ions. The PAD family members are thought to be involved in multiple sclerosis and rheumatoid arthritis pathophysiology, and they play a role in epidermis homeostasis (1). PAD6 is essential for formation of a novel oocyte-restricted fibrous structure, the cytoplasmic lattices (CPLs). PAD6/CPL superstructure plays a key role in regulating microtubule-mediated organelle positioning and movement (2).

References

- Vossenaar, E R. Et al: PAD, a growing family of citrullinating enzymes: genes, features and involvement in disease. *Bioessays*. 2003 Nov;25(11):1106-18.
- Kan, R. Et al: Regulation of mouse oocyte microtubule and organelle dynamics by PADI6 and the cytoplasmic lattices. *Dev Biol*. 2011 Feb 15;350(2):311-22.

Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **100 kDa**.

PAD6 Protein

Full length recombinant human protein expressed in Sf9 cells

Catalog #	P312-30FG
Lot #	T1065-3
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.