

PAD6 (Peptidyl arginine deiminase, type VI)

CATALOG NO.: PAD-21-494

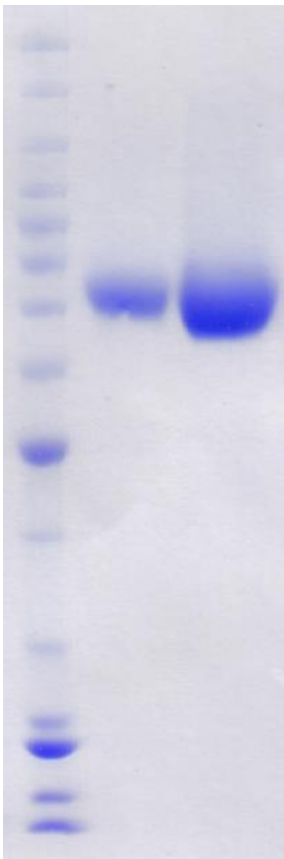
LOT NO.:

DESCRIPTION: Human recombinant PAD6 (residues 1-694; NCBI Reference Sequence NP_036519.2; MW = 80.1 kDa) expressed in *sf21* insect cells with an N-terminal Strep-tag. PAD6 (PADI6) catalyzes the hydrolytic deimination of specific arginine residues to citrulline and ammonia in a calcium-dependent manner.

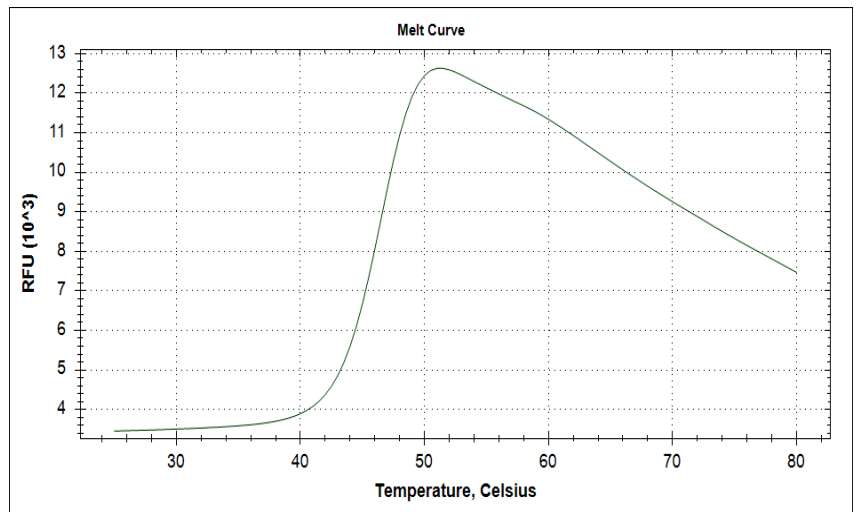
PURITY: >95% by SDS-PAGE

SUPPLIED AS: ___ µg/µL in 20 mM Tris, pH 8, 500 mM NaCl, 10% (v/v) glycerol, 1mM EDTA, 2mM DTT as determined by OD₂₈₀.

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue stained SDS-PAGE (4-12% acrylamide) of 4 and 10µg of RBC PAD6. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Differential Scanning Fluorimetry of RBC PAD6. Thermal denaturation of PAD6 is detected (CFX384 TMTouch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). The apo form of PAD6 displays a T_m of 46.5°C

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.