

Synthetic peptide: human PADI4

(Protein-arginine deiminase type-4)

Cat. No. C16000075

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide corresponding to the N-terminal region of human PADI4.

This peptide was used as immunogen to raise the Diagenode crude serum directed against PADI4 (Cat No. C15310075). The same peptide can be used to block this antibody in negative control experiments.

For more information about PADI4: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat No. C15310075).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: Aliquot upon receipt and store at -20°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Availability date: August 20, 2008

Last data sheet update: August 14, 2008