

PRODUCT NAME			
synthetic peptide: human PAPD5			
(PAP associated domain containing 5)			
Cat. No: sp-134-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide containing a sequence from the C-terminal part of human PAPD5
 This peptide was used as immunogen to raise the Diagenode crude serum directed against PAPD5 (cat# CS-134-100). The same peptide can be used to block this antibody in negative control experiments.
 For more information about PAPD5: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (cat# CS-134-100).

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: July 31, 2009

Last data sheet update: July 31, 2009

Lot #: 001/ Synthesis date: September 7, 2007/ aliquoting date: July 31, 2009