

Synthetic peptide: human PPARG

(peroxisome proliferator-activated receptor gamma)

Cat. No. C16000133

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide containing a sequence from the central part of human PPARG. This peptide was used as immunogen to raise the Diagenode crude serum directed against PPARG (Cat. No. C15410133). The same peptide can be used to block this antibody in negative control experiments. For more information about PPARG: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15410133).

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