

## Synthetic peptide: human IEF2C1/2

(eukaryotic translation initiation factor 2C, 1 and 2)

Cat. No. **C16000141**

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

**Description:** Synthetic peptide containing a sequence from the C-terminal part of human IEF2C1 and EIF2C2. This peptide was used as immunogen to raise the Diagenode crude serum directed against IEF2C1/2 (Cat. No. C15310141). The same peptide can be used to block this antibody in negative control experiments. For more information about IEF2C1/2: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15310141).

**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