

## Synthetic peptide: human ACTL6B

(actin-like 6B)

**Cat. No. C16000119**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide containing a sequence from the central part of human ACTL6B

This peptide was used as immunogen to raise the Diagenode crude serum directed against ACTL6B (Cat. No. C15310119). The same peptide can be used to block this antibody in negative control experiments.

For more information about ACTL6B: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15310119).

**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