

|                                    |                             |                               |                   |
|------------------------------------|-----------------------------|-------------------------------|-------------------|
| <b>PRODUCT NAME</b>                |                             |                               |                   |
| synthetic peptide:<br>human ACTL6B |                             |                               |                   |
| [actin-like 6B]                    |                             |                               |                   |
| <b>Cat. No:</b> sp-119-050         | <b>Format:</b> 50 µg/ 50 µl | <b>Concentration:</b> 1 µg/µl | <b>Lot #:</b> 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# CS-119-100). 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# CS-119-100).

**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: January 14, 2008/ aliquoting date: July 31, 2009