

<b>PRODUCT NAME</b> synthetic peptide: human p53			
<b>(Cellular tumor antigen p53)</b>			
<b>Cat. No:</b> sp-083-050	<b>Format:</b> 50 µg/ 50 µl	<b>Concentration:</b> 1 µg/µl	<b>Lot #:</b> 001

**Description:** Synthetic peptide corresponding to the C-terminal region of human p53.

This peptide was used as immunogen to raise the Diagenode ChIP grade antibody directed against p53 (cat# pAb-083-050). The same peptide can be used to block this antibody in negative control experiments.

For more information about p53: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (cat# pAb-083-050).

**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:** August 20, 2008

**Last data sheet update:** August 14, 2008

**Lot #:** 001/ Synthesis date: April 6, 2006/ aliquoting date: August 18, 2008