

# Progerin Monoclonal Antibody [13A4]

(Cat. No. A-0585)

## Background

Progerin is a malformed variant of lamin A, a protein vital in the nuclear lamina, a scaffold of proteins found inside the nuclear membrane of a cell. The genetic disorder Hutchinson-Gilford progeria syndrome (HGPS) is caused by a rare spontaneous mutation that leads to progerin being created in place of lamin A. Progerin is commonly generated by a silent point mutation (C1824T) in the lamin A gene (LMNA). This mutation triggers a cryptic splice site and gives rise to a form of lamin A with a deletion of 50 amino acids near the C-terminus. There is also evidence that small amounts of progerin is produced in normal cells and may be linked to the natural aging process.

## Description

Mouse monoclonal antibody to Progerin, Clone 13A4

## Formulation

Liquid. Neat cell culture supernatant containing 0.02% sodium azide.

## UniProt ID

P02545-6

## Immunogen

Synthetic KLH-coupled peptide corresponding to aa 604-611 (G604AQSPQNC611) at the C-terminus of human progerin.

## Specificity

Human

## Isotype

IgG1

## Purification

Serum

## Storage

Store at 4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

## Alternative Names

Delta50 Lamin A

## Application

ICC (1:10), IP (20 µl/assay), WB (1:1000)

## Ordering Information

### Products

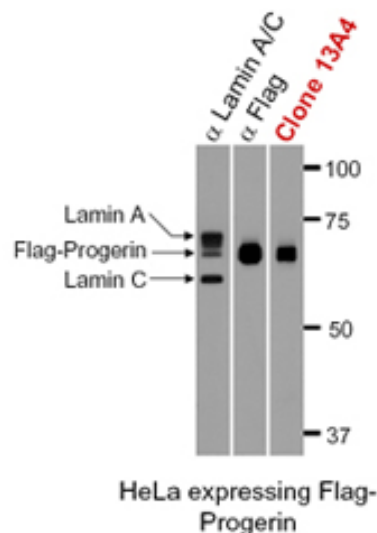
Progerin Monoclonal Antibody [13A4]

### Size

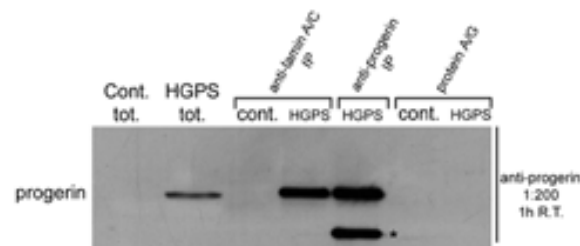
200 µl

### Cat. No.

A-0585-200



▲Western blot analysis (10% PAGE) of whole cell lysate of HeLa cells ectopically expressing FLAG-tagged human progerin. Lane 1: anti-Lamin A/C antibody (N18) (1:2000), Lane 2: anti-FLAG-tag antibody (M2) (1:5000), Lane 3: Progerin Monoclonal Antibody [13A4] (1:5)



▲Immunoprecipitation of progerin from lysates of Hutchinson-Gilford Progeria Syndrome (HGPS) cells. The immunoprecipitates were separated by SDS-PAGE, blotted and incubated with Progerin Monoclonal Antibody [13A4] (1:200). The asterisks labels the IgG band.

*This product is for research purposes only. Not intended for use in diagnostic procedures.*