

## AbFlex<sup>®</sup> Lamin A/C antibody (rAb)

**Catalog Nos:** 91189, 91190

**RRID:** AB\_2793795

**Isotype:** IgG2a

**Application(s):** IF, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 80 kDa

**Background:** AbFlex<sup>®</sup> antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports. AbFlex<sup>®</sup> Lamin A/C antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

Nuclear lamins are intermediate filament proteins that are the major structural component of the nuclear lamina on the inner surface of the nuclear envelope. Lamins A and Lamins C are splice variants of the Lamin A gene. Lamin A/C (CDCD1, LMN1, EMD2) expression is a hallmark of embryonic stem cell differentiation. In addition to adding structural integrity to the nucleus, lamins contribute to the makeup of the nuclear matrix. Lamins also help organize interphase chromatin through interactions with several chromatin proteins, including histones and Lap2, such that alteration in lamin organization results in disruption of DNA replication, transcription and RNA processing.

**Immunogen:** This Lamin A/C antibody was raised against a recombinant protein corresponding to amino acids 430-545 of human Lamin A.

**Buffer:** Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, 30% glycerol. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

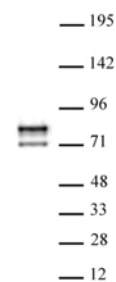
WB: 0.5 – 2 µg/ml

IF: 2 µg/ml

AbFlex<sup>®</sup> recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 39287.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

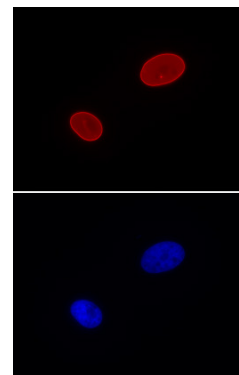
This product is for research use only and is not for use in diagnostic procedures.



AbFlex<sup>®</sup> Lamin A/C antibody (rAb) tested by Western blot.

10 µg of HeLa cell extract was probed with 1 µg/ml AbFlex Lamin A/C antibody.

MW: 80 kDa



AbFlex<sup>®</sup> Lamin A/C antibody (rAb) tested by Immunofluorescence

HeLa cells were stained with 2 µg/ml Lamin A/C antibody (upper panel) or DAPI (lower panel).