

Lamin A/C R482W antibody (mAb)

Catalog Nos: 39963, 39964

RRID: AB_2793411

Clone: 5H8-B4

Isotype: IgG1

Application(s): ICC, IF, IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 65 kDa

Background: 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 (laminopathy) results in disruption of DNA replication, transcription and RNA processing. The R482W mutation is involved in Dunnigan-type familial partial lipodystrophy (FPLD), a rare form of insulin resistance.

Immunogen: This Lamin A/C R482W antibody was raised against a recombinant protein corresponding to amino acids 430-545 of human Lamin A/C containing the R482W substitution.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91189.

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.