

## AIRE antibody (pAb)

**Catalog Nos:** 61139, 61140

**RRID:** AB\_2793523

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 60 kDa

**Background:** AIRE (Autoimmune regulator, APECED, Autoimmune polyendocrinopathy candidiasis ectodermal dystrophy) is a DNA-binding transcriptional regulator that promotes the expression of tissue-specific self-antigens in the thymus, which is important for self tolerance and the avoidance of autoimmune reactions. AIRE functions as a dimer or as a tetramer and interacts with nucleosomes, selectively interacting with histone H3 unmodified in the N-terminus through its first of two PHD domains. Defects in AIRE are a cause of autoimmune poly-endocrinopathy candidiasis ectodermal dystrophy (APECED) also known as autoimmune polyglandular syndrome type I (APS-1).

**Immunogen:** This AIRE antibody was raised against a peptide in the C-terminal region of human AIRE.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

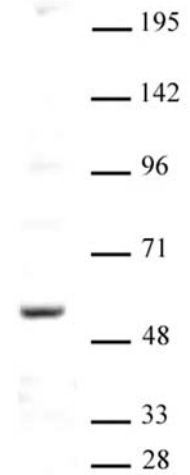
**Application Notes:**

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

**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.



**AIRE antibody (pAb) tested by Western blot.**

Detection of AIRE by Western blot. The analysis was performed using 30 µg Hep G2 whole-cell extract and AIRE (pAb) at a 1:500 dilution.