

PRODUCT INFORMATION

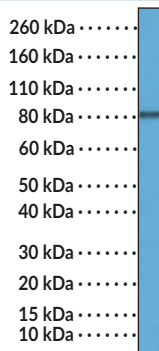
ALK/CD246 (C-Term) Rabbit Monoclonal Antibody (Clone RM361)

Item No. 32287

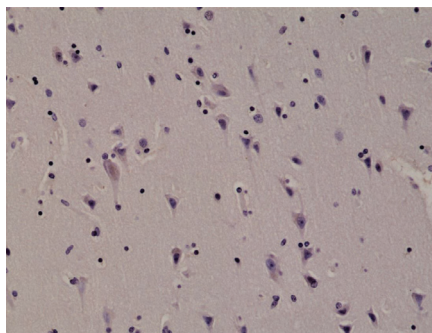
Overview and Properties

| | |
|----------------------------|---|
| Contents: | This vial contains 100 µl of protein A-affinity purified monoclonal antibody. |
| Synonyms: | ALK Tyrosine Kinase Receptor, Anaplastic Lymphoma Kinase, CD246, Cluster of Differentiation 246 |
| Immunogen: | Peptide from the C-terminal region of human ALK |
| Cross Reactivity: | (+) ALK, ALK fusion proteins |
| Species Reactivity: | (+) Human |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide |
| Clone: | RM361 |
| Host: | Rabbit |
| Isotype: | IgG |
| Applications: | Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

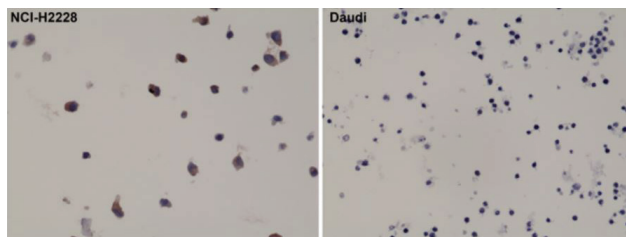
Images



WB of NSCLC cell line H2228 (cell lysate) expressing EML4-ALK variant 3 using ALK/CD246 (C-Term) Rabbit Monoclonal Antibody (Clone RM361) at a dilution of 1:2,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human brain tissue using ALK/CD246 (C-Term) Rabbit Monoclonal Antibody (Clone RM361) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded NCI-H2228 cells (expressing EML4-ALK variant 3) and Daudi cells (ALK negative) using ALK/CD246 (C-Term) Rabbit Monoclonal Antibody (Clone RM361).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/14/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Anaplastic lymphoma kinase (ALK), also known as CD246, is a receptor tyrosine kinase and member of the insulin receptor superfamily.¹ It is composed of an N-terminal extracellular domain that participates in cell-cell interactions, as well as transmembrane and juxtamembrane domains, and a C-terminal intracellular region that contains the protein kinase domain. Full-length ALK is expressed primarily in the embryonic nervous system, but the ALK kinase domain is expressed in a variety of cell types in the adult as a fusion protein.^{1,2} ALK is involved in cell proliferation and survival through downstream activation of the ERK1/2 and JAK/STAT signaling pathways, respectively.¹ The fusion protein nucleophosmin-ALK (NPM-ALK) has a constitutively active ALK kinase domain and is ectopically expressed in anaplastic large-cell lymphoma (ALCL), a type of non-Hodgkin lymphoma.³ A variety of ALK-based fusion proteins are associated with ALCL, non-small cell lung cancer (NSCLC), inflammatory myofibroblastic tumor (IMT), diffuse large B cell lymphoma (DLBCL), and other cancers.¹ Activating mutations in, and overexpression of, ALK have been associated with childhood neuroblastomas. Cayman's ALK/CD246 (C-Term) Rabbit Monoclonal Antibody (Clone RM361) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Roskoski, R., Jr. Anaplastic lymphoma kinase (ALK): Structure, oncogenic activation, and pharmacological inhibition. *Pharmacol. Res.* **68(1)**, 68-94 (2013).
2. Vernersson, E., Khoo, N.K.S., Henriksson, M.L., *et al.* Characterization of the expression of the ALK receptor tyrosine kinase in mice. *Gene Expr. Patterns* **6(5)**, 448-461 (2006).
3. Morris, S.W., Kirstein, M.N., Valentine, M.B., *et al.* Fusion of a kinase gene, ALK, to a nucleolar protein gene, NPM, in non-Hodgkin's lymphoma. *Science* **263(5151)**, 1281-1284 (1994).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM