

Anti-Cyclin E

Rabbit Polyclonal Antibody

Catalog # C82-363R

Lot # B3216-41

Cited Applications

ELISA, IP, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the 45kDa Cyclin E protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, Antiserum

Immunogen

The antibody was produced against synthesized peptide corresponding to a region near the carboxy terminus of human cyclin E conjugated to KLH.

Formulation

0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

Scientific Background

Anti-Cyclin-E antibody detects cyclin-E. Cyclin E is a member of the cyclin family. Cyclin E binds to the G1 phase Cdk2, which is required for the transition from G1 to S phase of the cell cycle that determines cell division. The Cyclin E/CDK2 complex phosphorylates p27Kip1, tagging it for degradation, thus promoting expression of Cyclin A, allowing progression to S phase. Anti-cyclin-e Antibody is ideal for investigators involved in Cell Signaling, cell biology and Signal Transduction research.

References:

1. Koff, A., et al.: Human cyclin E, a new cyclin that interacts with two members of the CDC2 gene family. *Cell*. 1991: 66; 1217-1228.
2. Koff, A., et al.: Formation and activation of a cyclin E-cdk2 complex during the G1 phase of the human cell cycle. *Science*. 1992: 257; 1689-1694.
3. Claudio, P.P., et al.: Functional analysis of pRb/p130 intereaction with cyclins. *Cancer Res*. 1996: 56; 2003-2008.

Anti-Cyclin E

Rabbit Polyclonal Antibody

Catalog #	C82-363R
Lot #	B3216-41
Purification	Delipidation and defibrination
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.