

Anti-DYKDDDDK

Rabbit Polyclonal Antibody

Catalog # F51-63R

Lot # R027-7

Cited Applications

- Western blot (1:5000 dilution)

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

Specificity

Recognizes the DYKDDDDK tag

Cross Reactivity

- All proteins tagged with DYKDDDDK epitope

Host

Rabbit

Immunogen

DYKDDDDK peptide conjugated to KLH

Formulation

TBS, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

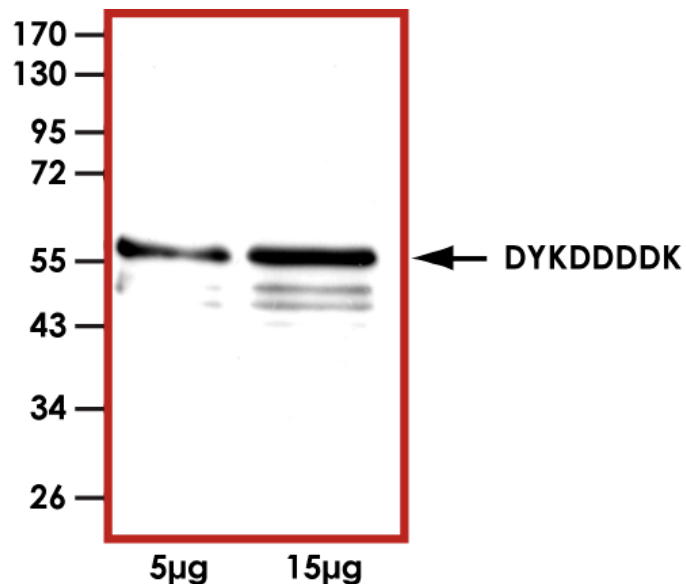
Scientific Background

The DYKDDDDK octapeptide is small, highly charged, and hydrophilic, making it useful across a wide variety of applications [1]. It is commonly employed in structural and functional studies where it is necessary to preserve the native conformation and and/or activity of the protein [2].

References

1. Jarvik, JW. and Telmer, CA: Epitope tagging, Annual Review of Genetics, 1998; 32, 601-618.
2. Larson DM. et al: Functional expression and biochemical characterization of an epitope-tagged connexin37. Mol Cell Biol Res Commun 3: 2000; 115-121.

Sample Data



Representative western blot with Anti-DYKDDDDK (1:5000) using 5 µg, and 15 µg of a positive cell lysate transfected with a DYKDDDDK tagged fusion protein.

Anti-DYKDDDDK

Rabbit Polyclonal Antibody

Catalog Number

F51-63R

Specific Lot Number

R027-7

Purification
Concentration

Affinity chromatography
1.0 µg/µL

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.