

Dehydrin (plant) polyclonal antibody

Dehydrins are a family of proteins that become abundant during dessication in seedlings and embryos of cereal crop plants including barley, corn, wheat, and rice. Dehydrin proteins have a highly conserved lysine-rich block (KIKEKLPG) found near the carboxy terminus which appears to be characteristic of, and unique to Dehydrins.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-PLA-100-D	50µg
ADI-PLA-100-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Dehydrin (plant), pAb: Lane 1: MW marker, Lane 2: Corn.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application Notes Detects a band of ~65kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID Y00842 (rice dehydrin family member Rab21)

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of dehydrin. The sequence is completely conserved in barley, rice, maize and wheat.

Purity Detail Protein A affinity purified.

Recommendation Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Plant

UniProt ID A2ZDX9 (rice dehydrin family member Rab21)

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com