

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 Dilutions/Conditions	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](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

Last modified: May 29, 2024