

DGPx2 (Deamidated Gliadin Peptide – fusion protein)

DGPx2 is a fusion protein of two different deamidated gliadin peptides, fused with a carrier protein to be used as antigen for the detection of antibodies specific for deamidated gliadin.

The recombinant antigen is meant for solid (ELISA and immunoblot) and fluid phase diagnostic assays. The protein is bound by human type IgA and IgG (auto) antibodies.

Ordering Information [Order Online »](#)

NZY-G006-001	1mg
NZY-G006-250	250µg

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Handling	After reconstitution, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White lyophilized solid.
Formulation	Lyophilized from a solution of ~50mM sodium dihydrogen phosphate, pH 6.8.
MW	~31kDa
Purity	≥95% (SDS-PAGE and Coomassie staining)
Reconstitution	Add at least the volume of water the protein is lyophilized with (see Certificate of Analysis) to the vial of lyophilized powder. Rotate vial gently until solid dissolves. After reconstitution the solution should be stored frozen in working aliquots.
Source	Produced in <i>E. coli</i> .
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



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