

PDQ Protease Assay™

The PDQ Protease Assay is a colorimetric assay used to detect protease activity. The proprietary substrate is susceptible to a wide range of proteases, e.g., collagenase, proteinase K, papain, pepsin, bromelin, ficin, trypsin and chymotrypsin. This substrate is a cross-linked matrix containing protein and a dye-protein conjugate. The matrix is supplied in ready-to-use vials. Test samples containing putative protease activity are aliquoted into the vials which are then incubated at an appropriate temperature. Protease activity is detected spectrophotometrically with increasing optical density proportional to increasing enzyme activity.

Ordering Information

[Order Online »](#)

AES-0201	1Kit
----------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents 48 vials containing 200µl of pre-made matrix

Technical Info / Product Notes Buffers that can be used to prepare test samples may range in pH from 4 to 10.

Materials not provided but required:

– 0.2N NaOH

– Spectrophotometer for reading absorbance at 450 nm

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