

Sample Deproteinizing Kit (PCA)

To remove proteins from various sample types.

It is a fundamental requirement in biochemistry and related fields to be able to remove those constituents of the sample under investigation which will interfere with analysis in any of several ways, both physical and chemical. When the analyte in question is a small molecule, the presence of macromolecules such as proteins, nucleic acids and carbohydrates can seriously impair analysis. Proteins are usually removed by precipitation using one of several chemicals including acetone, perchloric acid (PCA), metaphosphoric acid, sulfosalicylic acid or trichloroacetic acid. PCA is a strong acid commonly used for protein precipitation and removal in biological samples. AkrivisBio's Deproteinizing Sample Preparation Kit utilizes a PCA precipitation method, providing a unique tool to prepare samples for the analysis of a variety of small molecules such as lipids, metabolites etc.

Ordering Information

[Order Online »](#)

AKR-PI-0103	200preparations
-------------	-----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Species independent

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