

Product datasheet

FASP Protein Digestion Kit ab270519

18 References

Overview

| | |
|-------------------------|---|
| Product name | FASP Protein Digestion Kit |
| Product overview | <p>FASP Protein Digestion Kit (ab270519) is for researchers who wish to solubilize whole or fractionated protein samples in SDS, digest the protein with trypsin, and analyse the resulting peptides by mass spectrometry. The FASP Protein Digestion Kit provides the necessary columns and buffers to carry out Universal Sample preparation as developed by Wisniewski, Zoubman, Nagaraj and Mann (2009). Ab270519 is compatible with a comprehensive range of biological sample types.</p> <p>This product is manufactured by Expedeon, an Abcam company and it was previously called FASP Protein Digestion Kit. Expedeon product code 44250 is the same as the 8-pack size of this product.</p> |

Properties

| | |
|-----------------------------|---|
| Storage instructions | Store at room temperature. Please refer to protocols. |
|-----------------------------|---|

| Components | 8 packs |
|--|------------|
| ab274283 - 0.5 M Sodium Chloride Solution | 1 x 1ml |
| ab274284 - 50 mM Ammonium Bicarbonate Solution | 1 x 20ml |
| ab280352 - Collection tube | 16 units |
| ab274286 - Iodoacetamide | 8 x 9800µg |
| ab274288 - Spin Filter | 8 units |
| ab274285 - Tris Hydrochloride Solution | 1 x 20ml |
| ab274287 - Urea | 8 x 750mg |

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit [**https://www.abcam.com/abpromise**](https://www.abcam.com/abpromise) or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors