

Product Information

Fatty Acid Methyl Ester Purification Kit

Manufactured by Nacalai Tesque, Inc.

Catalog Number **MAK225**
Store at Room Temperature

TECHNICAL BULLETIN

Product Description

The Fatty Acid Methyl Ester Purification Kit is used to further purify samples prepared with the Fatty Acid Methylation Kit (Catalog Number MAK224). Methylated fatty acid samples are required to be purified with this kit before the analysis by a capillary column to prevent premature column deterioration caused by the reaction impurities.

Components

The kit is sufficient for 50 sample purifications.

Conditioning Solution Catalog Number MAK225A	200 mL
Washing Solution Catalog Number MAK225B	200 mL
Eluting Solution Catalog Number MAK224C	200 mL
SPE Cartridge Column Catalog Number MAK225D	50 ea

Reagents and Equipment Required but Not Provided.

- Fatty Acid Methylation Kit, Catalog Number MAK224

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

The kit is shipped at ambient temperature. Store all components at room temperature upon receiving.

Procedure

All samples and standards should be run in duplicate.

All steps are done under gravity fall condition.

- Add 3 mL of Conditioning Solution to a SPE Cartridge Column.
- Add the methylated sample from the Fatty Acid Methylation Kit (Catalog Number MAK224) to the SPE Cartridge Column.
- Add 3 mL of Washing Solution to the SPE Cartridge Column.
- Add 3 mL of Eluting Solution and collect eluted liquid containing fatty acid methyl esters.

SJ,MAM 03/16-1