

# PRODUCT INFORMATION



## SPE Cartridges (C18) (1 ml)

Item No. 400077

### Overview

---

**Synonyms:** C18 Column, Solid Phase Extraction Cartridges  
**Storage:** Room temperature  
**Stability:** As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly

### Laboratory Procedures

---

This product may be used for purification of various lipid mediators such as prostaglandins and hydroxylated fatty acids from various sample types prior to analysis by LC-MS/MS or ELISA. References are provided below for applicable C18 sample purification methods prior to LC-MS analysis.<sup>1-4</sup> For sample purification prior to analysis by ELISA, we recommend consulting the appropriate assay kit booklet for instructions.

### Precaution

---

This product may not perform as described if used with other assay reagents or protocols.

### References

---

1. Maddipati, K.R. and Zhou, S.-L. Stability and analysis of eicosanoids and docosanoids in tissue culture media. *Prostaglandins and other Lipid Mediators* **94**, 59-72 (2011).
2. Maddipati, K.R., Romero, R., Chaiworapongsa, T., et al. Eicosanomic profiling reveals dominance of the epoxygenase pathway in human amniotic fluid at term in spontaneous labor. *FASEB J.* **28(11)**, 4835-4846 (2014).
3. Powell, W.S. Rapid extraction of arachidonic acid metabolites from biological samples using octadecylsilyl silica, Chapter 58, in *Methods in Enzymology*. Academic Press, New York, 467-477 (1982).
4. Song, W.-L., Wang, M., Ricciotti, E., et al. Tetranor PGDM, an abundant urinary metabolite reflects biosynthesis of prostaglandin D<sub>2</sub> in mice and humans. *J. Biol. Chem.* **283(2)**, 1179-1188 (2008).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/13/2019

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM