

Agilent Captiva EMR—Lipid Method Guide

General instructions for 96-well plate and 1 mL cartridge formats:

Agilent Captiva EMR—Lipid cartridges and 96-well plates allow streamlined in-well protein precipitation, filtration, and cleanup for lipid-containing samples. The improved filter design gives easy flow with vacuum or positive pressure, and allows for protein precipitation without clogging during elution. The novel EMR—Lipid sorbent chemistry provides highly selective and efficient lipid/matrix removal without impacting analyte recovery. Effective matrix removal assures minimal ion suppression or enhancement on target analytes, which significantly improves method reliability and ruggedness.

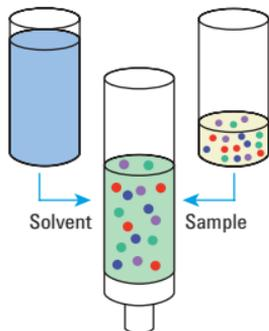
The 96-well plates are ideal for high-throughput workflows, while 1 mL cartridges can accommodate small batch needs.



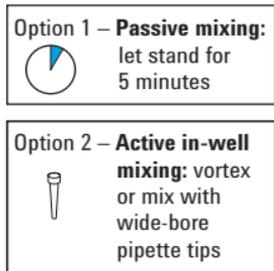
Agilent Technologies

Operating instructions and tips for Agilent Captiva EMR—Lipid 96-Well Plate and 1 mL Cartridge Products

1. Add biological sample and crash solvent



2. Mix to precipitate protein



3. Filter



It is highly recommended to add sample first and then crash solvent, to achieve better sample homogeneity during sample and solvent addition.

● Salts ● Proteins ● Lipids ● Analyte



User Tips

Protein precipitation workflow

Product configuration	96-well plate or 1 mL cartridge
Sample size	Between 20–200 μ L
Sample treatment	Crash solvent ratio: between 3:1 and 5:1 ACN with 1 % formic acid to sample. If total volume is less than 500 μ L, add additional 4:1 ACN:H ₂ O to reach a minimum volume of 500 μ L. ACN is preferable to MeOH to maximize protein precipitation and avoid gelation.
Recommended sample addition order	1) Sample 2) Crash solvent
Mixing	Option 1: Passive mixing. Let stand for 5 minutes to allow for complete protein precipitation to occur. Option 2: Active in-well mixing. For 96-well plates containing >500 μ L of total volume and 1 mL cartridges regardless of volume, pipette mixing using wide bore pipette tips (perform 3 to 5 aspiration/dispense cycles) is recommended. For 96-well plates with <500 μ L of total volume, cap and vortex at 1,350 rpm. Option 3: Protein precipitation and mixing can be performed in a separate tube, centrifuged, and subsequently transferred to the Captiva EMR—Lipid well/cartridge.
Pass-through filtration and cleanup	Vacuum between 2–5 in Hg initiates flow. Positive pressure (3–4 psi) is also acceptable. For optimal lipid removal, a controlled flow rate of one drop every 3–5 seconds is highly recommended. After elution, apply higher vacuum or positive pressure for 1–2 minutes to ensure maximum sample recovery. Flow rate is dependent on sample type, crash solvent, and mixing. An alternative approach to vacuum and positive pressure is centrifugation. For 96-well plates, 500–800 rpm for a minimum of 10 minutes is recommended, followed by 1–2 minutes at 2,500–3,000 rpm to drain the cartridge. Centrifugation speed and time are dependent on the sample volume and matrix.

For more detailed protocol recommendations, please refer to Agilent Application Note 5991-9222EN.

Agilent Captiva EMR—Lipid ordering information

Part number	Description	Quantity
5190-1000	Agilent Captiva EMR—Lipid 96-well plate	1 plate
5190-1001	Agilent Captiva EMR—Lipid 96-well plate	5 plates
5190-1002	Agilent Captiva EMR—Lipid 1 mL cartridge	100/pk

Learn more
agilent.com/chem/samplepreparation

Buy online
www.agilent.com/chem/store

Find an Agilent office or authorized distributor
www.agilent.com/chem/contactus

USA and Canada
1-800-227-9770
agilent_inquiries@agilent.com

Europe
info_agilent@agilent.com

Asia Pacific
inquiry_lsca@agilent.com

India
india_lsca_marketing@agilent.com

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2017
Printed in the USA, August 23, 2018
5991-7930EN



Agilent Technologies