

# MARATHON FILAMENT FOR PERKINELMER GC/MS SYSTEMS

**Marathon Filament**  
a revolutionary technology developed exclusively  
by PerkinElmer, delivering long life even under  
the most difficult chromatography conditions.

After lengthy performance testing and filament research, the Marathon Filament has been engineered to provide exceptional long life and withstand difficult chromatography conditions. It has high resistance to demanding injections such as headspace or purge and trap and has stood up to challenging applications such as flame-retardant analysis.

The Marathon Filament is a direct replacement of the previous rhenium filament – no system changes required.

At PerkinElmer, we understand your challenges and are committed to providing the best solutions to make your life easier – the Marathon Filament is part of our ongoing efforts to deliver the latest technology for our PerkinElmer GC/MS platform.

Part No.	Description
N6470012	Marathon Filament for PerkinElmer GC/MS Systems

To view the webcast and to order your replacement filament, visit [www.perkinelmer.com/MarathonFilament](http://www.perkinelmer.com/MarathonFilament)

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)

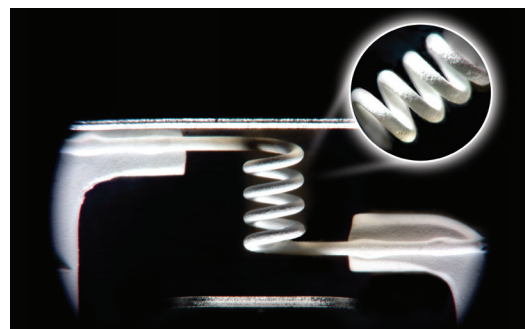
  
**PerkinElmer**<sup>®</sup>  
For the Better

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2015, PerkinElmer, Inc. All rights reserved. PerkinElmer<sup>®</sup> is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

012291\_01

PKI



## Features and Benefits

- Long life even under the most difficult chromatography conditions
- Unique white surface engineered for maximum durability and optimum performance
- Now included with all new Clarus<sup>®</sup> GC/MS systems (560 and 600 series)
- Backward compatible with all units using rhenium filaments
- Works with both electron and chemical ionization sources