

Instruction Manual Supplement

说明书补充

Manual Title: Model 5030 Synchronized Hybrid Ambient Real-time Particulate Monitor

说明书名称: 型号 5030 同步混合环境空气实时颗粒物监测仪

Part No: 102017-00

说明书编号: 102017-00

Issue date: 8Jan2013

版本日期: 2013 年 1 月 8 日

Revision date: 5Feb2018

修改日期: 2018 年 2 月 5 日

This supplement contains information necessary to ensure the accuracy of the above manual. This manual is distributed as an electronic manual on <https://www.thermofisher.com/> For more information, Please consult your local sales representative for details.

本说明书补充载有必要的信息以确保上述说明书的准确性。本说明书以电子版本在 <https://www.thermofisher.com/> 发行。有关详细信息，请咨询当地销售代表。

Change #1

On page 17, add the following under the 2nd paragraph:

在第 7 页第二自然段后增加以下说明：

The %RH threshold setting range of 30% to 60% is suggested. %RH threshold setting and minimum heating power should be properly set with consideration of the local meteorological conditions and particulate matter composition. This setting should be compatible with local regulatory agency considerations and should be set based on comparison with gravimetric sampling method results at the same location as the instruments being operated. Always refer to local regulations prior to setting this value.”

RH% 阈值设置范围建议在 30%-60%。在考虑当地气象条件和颗粒物组成的情况下，用户应适当设置 RH% 阈值以及最小加热功率。此设置应与当地监管机构的意见相一致，并应根据与仪器运行位置相同的手工监测方法（重量法）的分析结果进行设置。在参数设置时，请始终参考当地法规。

Change #2

On page 35, add the following under Table 2-1. EU and US monitoring protocols:

在第 35 页表 2-1 EU 和 US 监测协议下增加以下说明：

The protocols below are only example of comparison between EU and US. For China and other countries, the %RH threshold setting range of 30% to 60% is suggested. The value should be set based on consideration of the local meteorological conditions and particulate matter composition. The setting should be based on a comparison with gravimetric sampling method results at the same location as the instruments being operated.” Always refer to local regulatory agencies for guidance prior to setting this value.

下述设置方案仅是欧盟和美国的应用案例。对于中国和其他国家，RH% 阈值设置范围建议在 30%-60%。该设置值应考虑到当地气象条件和颗粒物组成并应根据与仪器运行位置相同的手工监测方法（重量法）的分析结果进行设置。在参数设置时，请始终参考当地法规。