

# PDM3700 Battery Charger Update

Thermo Fisher Scientific has released a new version of the charger (PN 57-010925-3700) starting with Serial Number 2006489. Charging using the new version is performed in the same way as with the original. This technical note describes the charging and front panel LED function of the updated charger for use with the PDM3700, some of which may slightly differ from the original charger:

## LED Function

- **LED Off:** Indicates that the power to the charger is off or no PDM is connected to the charger.
- **LED Solid Red:** Battery is being charged and the elapsed charge time is less than 7 hours.
- **LED Flashing Green (Fast, 1 Hz):** The battery is being charged and the charge time has exceeded 7 hours. The battery charge level is at least 80%.
- **LED Flashing Green (Slow, 0.5 Hz):** The battery charge time has reached 14 hours. The integrated circuit within the charger has not indicated that the battery is fully charged, but the battery voltage and charge time indicate that the battery is essentially fully charged.
- **LED Solid Green:** This will occur after a charge time of less than 14 hours and indicates that the battery pack is fully charged.
- **LED Flashing Red:** There are three conditions that trigger a Flashing Red condition:
  1. Battery voltage does not reach a minimum threshold voltage level within 60 minutes.
  2. Battery voltage is below a threshold limit after 7 hours will not fully charge in the maximum charge time, indicating a fault with the battery pack or charging cycle.
  3. The charger detects a fault condition, such as, low voltage, over voltage, or high battery temperature. Contact Thermo Fisher Scientific Technical support for assistance.

## Notes:

- At the end of a charge cycle with a Solid Green LED, the LED may briefly turn to solid red or flashing green conditions. This is normal behavior and does not indicate a fault with the battery pack or charger.
- If the charger is turned off and back on, or if the charger cable is removed and reconnected, the charging cycle could restart as outlined above. Typically, 30 – 40 mins are required to allow the circuit to detect the current charge state.
- If the PDM3700 is not being recognized, reconnecting the charger cable will correct this condition.

27 Forge Parkway  
Franklin, MA 02038  
Ph: (508) 520-0430  
Toll Free: (866) 282-0430  
orders.aqi@thermofisher.com

Find out more at [thermofisher.com](https://www.thermofisher.com)

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified

**ThermoFisher**  
SCIENTIFIC