

NanoDrop Supplemental Warranty

Thermo Fisher Scientific warrants that each NanoDrop™ spectrometer we sell is free from defects in labor and materials and shall conform to its product specifications as defined in the product user documentation. If the product does not function as warranted during the warranty period, we will repair or replace it without charge. If in our judgment we are unable to do so, you may return it to us and we will refund your money.

This warranty replaces all other warranties, expressed or implied, including the implied warranties of merchantability and fitness for a particular purpose and any other obligations or liabilities on our part whether in contract, warranty, negligence or otherwise. We shall not be liable for and disclaim all consequential, incidental and contingent damages.

Warranty Period

The warranty period is 24 months. If you install the product, the warranty begins on date of invoice. If we install the product, the warranty period begins on the date of installation but will begin no later than 30 days from the date of the invoice.

The warranty period for products sold outside the U.S.A. and Canada is 24 months from the date of installation or 26 months from the date of shipment, whichever is less.

Limit of Warranty

Misuse, accident, modification, unsuitable physical or operating environment, improper maintenance, or damage caused by a product for which we are not responsible will void the warranty.

Items Not Covered by Warranty

We do not warrant uninterrupted or error-free operation of the product.

We provide certain non-Thermo Fisher Scientific products on an “as is” basis. Non-Thermo Fisher Scientific manufacturers or suppliers may provide their own warranties to you.

Certain components may have separate warranty periods as stated in the product user documentation. Consumables are not covered by warranty.

A separate software warranty is provided with the software end user license agreement.