

**Date:** January 30, 2020

**Subject:** Notification - Obsolescence of Nuclear Level and Density Gauges

At Thermo Fisher Scientific, we continually evaluate our global operations and product portfolios to identify opportunities to improve efficiency and effectiveness in meeting the needs of our global customers.

After a careful and thorough review of our Environmental and Process Monitoring business, we have decided to discontinue our Nuclear Level and Density product line, effective January 31, 2020.

The products impacted by this decision are listed below:

<b>Product Description</b>	<b>Product Model</b>	<b>Replacement</b>
<b>Gamma density gauge</b>	<b>DensityPRO (MS2011DU, MS2011DUR)</b>	<b>None</b>
<b>Gamma level gauge</b>	<b>LevelPRO (MS2011LU1, MS2011LU1R)</b>	<b>None</b>
<b>NAI density gauge</b>	<b>DensityPRO NAI (NS2011I, MS2011R)</b>	<b>None</b>
<b>3680 Smart density transmitter</b>	<b>3680, CutPRO, ConsistencyPRO</b>	<b>None</b>
<b>InterfacePRO for level, density, and interface</b>	<b>InterfacePRO (MS2011IP)</b>	<b>None</b>
<b>DensityPRO integrated gamma system</b>	<b>DensityPRO-C</b>	<b>None</b>
<b>SGD-O density gauge for fracturing</b>	<b>SGDO Density Gauge</b>	<b>None</b>
<b>KRILPRO neutron backscatter level/interface device</b>	<b>KRILPRO (7200A)</b>	<b>None</b>
<b>MOLA-Moisture Online Analyzer</b>	<b>MOLA/MOLA-LS (7200A)</b>	<b>None</b>
<b>PNF Point level gauge</b>	<b>PNF Point Level Switch</b>	<b>None</b>

Thermo Fisher will continue to provide service and support for Nuclear Level and Density products, subject to availability, through the end of December 2024.\*

The duration is based on parts availability and begins with the last date of manufacture or declared product obsolescence date, whichever is earlier. Thermo Fisher Scientific reserves the right to change the level or length of support without notice.

You are a valued and important customer and we are committed to supporting you through this transition. We understand you may have questions or concerns regarding this decision. Please be advised that your Thermo Fisher partner can support you through this change. Please reach out to them for additional support.

Sincerely,

**Ben Fuentes**

Product Line Manager

Nuclear Level and Density

Thermo Fisher Scientific

**\*It is possible that certain consumables and spare parts could become unavailable prior to the end of the defined period due to supplier availability or regulatory restrictions.**