

COMPREHENSIVE SERVICE SOLUTIONS

KRILPRO Upgrade Kit

Upgrade Today

Get **10% off***
your total order

- 10% discount applies to KRILPRO upgrade kit plus Field Service, Contracts and/or Spare Parts
- Now thru 12/31/2020
- Must mention discount code: 2020 **Service KRILPRO Upgrade**
- Certain restrictions apply, see your Sales Rep for details

UPGRADE



The discontinuation of previous generation instruments is part of our goal to continuously offer you our latest technology and the best solutions to meet your measurement needs. The legacy KRIL is obsolete and no longer supported. Upgrade today to ensure long term uptime with supported maintenance and spare parts inventory.

Powered by an extremely small source, the Thermo Scientific™ KRILPRO fills the need for an ALARA-compliant, yet highly precise, level/interface detection system. Rugged and reliable, it features neutron backscatter technology and is engineered to endure harsh environments to help optimize delayed coking operations and improve profitability.

Features

- KRILPRO upgrade allows you to re-use existing Am-Be source
- Measurement through thick-walled vessels of any diameter
- Non-contacting measurement unaffected by temperature, pressure, viscosity, corrosives or abrasives
- Continuous self-diagnostics provide instant check of system integrity
- Requires minimal maintenance and ensures no unnecessary downtime
- Precision of less than 0.5% of span

USA

27 Forge Parkway
Franklin, MA 02038
Ph: (713) 272-0404
Fax: (713) 272-2273
orders.process.us@thermofisher.com

Europe

Ion Path, Road Three,
Winsford, Cheshire CW73GA UK
Ph: +44 1606 548700
Fax: +44 1606 548711
sales.epm.uk@thermofisher.com

*Can not be combined with other promotions or contracted discounts. Not valid on previous purchases. Must purchase the KRILPRO upgrade kit at the same time as field service, contracts and/or spare parts. Discount only applicable to sales made with a binding purchase order received on or before 12/31/2020. Customers must accept delivery by 12/31/2020.

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