



## Certificate of Registration of Quality Management System to ISO 9001:2015

The National Standards Authority of Ireland Inc. certifies that:

**Revvity**  
**2650 Crescent Drive**  
**Suite 200**  
**Lafayette, CO 80026**  
**USA**

has been assessed and deemed to comply with the  
requirements of the above standard in respect of the scope of  
operations given below:

---

**Manufacture, kit, and distribution of molecular  
biology reagents and kits for research use only.**

Geraldine Larkin  
Chief Executive Officer

Lisa Greenleaf  
Technical Operations Officer

---

Registration Number: 19.8127  
Certification Granted: August 24, 2017  
Effective Date: February 03, 2025  
Expiry Date: February 02, 2028



---

National Standards Authority of Ireland Inc., 20 Trafalgar Square, Nashua, New Hampshire, NH 03063, USA T +1 603 882 4412