



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

The National Standards Authority of Ireland Inc. certifies that:

Revvity Health Sciences Inc.
549 Albany Street
Boston, MA 02118
USA

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

**Process Development and Manufacture of Radio
Labeled Specialty Chemicals, Non-labeled
Analogues, Radionuclides, Diagnostic and Research
Detection Systems (kits) and Reagents, Molecular
Biology Kits and Reagents, and Custom
Membranes, Kits and Reagents**

Geraldine Larkin
Chief Executive Officer

Lisa Greenleaf
Technical Operations Officer

Registration Number: 19.4411
Certification Granted: February 19, 2009
Effective Date: November 18, 2024
Expiry Date: November 17, 2027

