

Licensing Number **12759**

Licence Holder **Revvity (UK) Ltd**

Licensed Premises **Revvity (UK) Ltd**
Suite 23LG05-23, Block 23
Alderley Park
Congleton Road
Nether Alderley
Macclesfield
SK10 4TG

Designated Individual **Mr Umar Farooq**

This licence is granted under Section 16 (2) (e) (ii) of the Human Tissue Act 2004.

This licence authorises the storage of relevant material which has come from a human body for use for the following scheduled purposes:

- Determining the cause of death
- Establishing after a person's death the efficacy of any drug or other treatment administered to him
- Obtaining scientific or medical information about a living or deceased person which may be relevant to any other person (including a future person)
- Public display
- Research in connection with disorders, or the functioning, of the human body
- Clinical audit
- Education or training relating to human health
- Performance assessment
- Public health monitoring
- Quality assurance

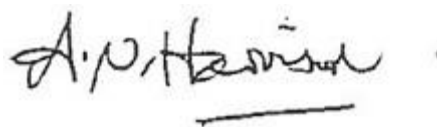
The licensed activity should be carried on only at the licensed premises specified above, and under the supervision of the Designated Individual.

This licence is subject to the conditions set out in the Annexes accompanying this licence as may be subsequently varied pursuant to an application under paragraph 8 of Schedule 3 to the Human Tissue Act 2004.

This licence is valid from the date specified below and will remain in force until revoked.



.....
Lynne Berry CBE
Chair



.....
Nicolette Harrison
Director of Regulation

Valid from: 24 June 2025

Annex A – Additional conditions

**Revvity (UK) Ltd
Suite 23LG05-23, Block 23
Alderley Park
Congleton Road
Nether Alderley
Macclesfield
SK10 4TG**

There are no additional conditions applied to this licence.