



# NSAI

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

The National Standards Authority of Ireland Inc. certifies that:

**Revvity Health Sciences Private Limited**  
**Block No. 702 – 704, 7th Floor**  
**TICEL Bio Park (Phase II)**  
**No.5, CSIR Road, Taramani**  
**Chennai, 600113**  
**Tamil Nadu, India**

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

---

### **Screening and Confirmatory Tests that Relate to Indicators for Human Genetic Diseases Focused on Fetal, Maternal and Newborn Domain**

**Additional sites covered under this multi-site certification are listed on the Annex (File No. MD19.4828)**

Geraldine Larkin Chief  
Executive Officer

Lisa Greenleaf  
Technical Operations Officer

---

Registration Number: 19.4828  
Certification Granted: June 23, 2015  
Effective Date: June 23, 2024  
Expiry Date: June 22, 2027





# NSAI

## **Annex to Certificate Number: 19.4828**

### **Scope of Registration:**

**Screening and Confirmatory Tests that Relate to Indicators for Human Genetic Diseases Focused on Fetal, Maternal and Newborn Domain**

#### **Activity**

#### **Location**

Central Management Group,  
Prenatal and Neonatal Screening,  
Confirmatory Tests

Revvity Health Sciences Private Limited  
Block No. 702 – 704, 7th Floor  
TICEL Bio Park (Phase II)  
No.5, CSIR Road, Taramani  
Chennai, 600113  
Tamil Nadu, India  
File No.: 19.4828

Prenatal Screening

Revvity Health Sciences Private Limited  
3-5-874/1, Opp. Toyota Showroom,  
Hyderguda, Hyderabad, Telangana  
500029  
India  
File No.: 19.4828/B

**Verified by:  
Technical Operations Officer**