



NSAI

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

The National Standards Authority of Ireland Inc. certifies that:

Revivity 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: October 25, 2023
Expiry Date: June 22, 2024





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**