



National Accreditation Board for
Testing and Calibration Laboratories

CERTIFICATE OF ACCREDITATION

REVVITY HEALTH SCIENCES PRIVATE LIMITED

has been assessed and accredited in accordance with the standard

ISO 15189:2012

**"Medical laboratories - Requirements for quality and
competence"**

for its facilities at

BLOCK NO. 702-704, 7TH FLOOR, TICEL BIO PARK, PHASE II, CSIR RD, THARAMANI, CHENNAI, TAMIL
NADU, INDIA

in the field of

Medical Testing

Certificate Number: MC-6167

Issue Date: 04/12/2023

Valid Until:

22/11/2025

This certificate remains valid for the Scope of Accreditation as specified in the annexure subject to continued
satisfactory compliance to the above standard & the relevant requirements of NABL.
(To see the scope of accreditation of this laboratory, you may also visit NABL website www.nabl-india.org)

Name of Legal Entity: REVVITY HEALTH SCIENCES PRIVATE LIMITED

Signed for and on behalf of NABL



N. Venkateswaran
Chief Executive Officer



National Accreditation Board for Testing and Calibration Laboratories

SCOPE OF ACCREDITATION

Laboratory Name :

REVVITY HEALTH SCIENCES PRIVATE LIMITED, BLOCK NO. 702-704, 7TH FLOOR, TICEL BIO PARK, PHASE II, CSIR RD, THARAMANI, CHENNAI, TAMIL NADU, INDIA

Accreditation Standard

ISO 15189:2012

Certificate Number

MC-6167

Page No

1 of 3

Validity

04/12/2023 to 22/11/2025

Last Amended on

-

S.No	Discipline / Group	Materials or Products tested	Component, parameter or characteristic tested / Specific Test Performed / Tests or type of tests performed	Test Method Specification against which tests are performed and / or the techniques / equipment used
Permanent Facility				
1	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Biotinidase	Time Resolved Fluoroimmunoassay
2	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Free Beta - Human Chorionic Gonadotropin (Free β -hCG)	Time Resolved Fluoroimmunoassay
3	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Neonatal 17-a- Hydroxy-progesterone (17OHP)	Time Resolved Fluoroimmunoassay
4	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Neonatal Glucose - 6 - Phosphate Dehydrogenase (G6PD)	Oxidation of Glucose - 6 - Phosphate
5	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Neonatal Immunoreactive Trypsin(ogen) (IRT)	Time Resolved Fluoroimmunoassay
6	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Neonatal Thyroid Stimulating Hormone (hTSH)	Time Resolved Fluoroimmunoassay
7	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Neonatal Total Galactose (TGal)	Fluorescence, Galactose Oxidase Method
8	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Phenylalanine	Ninhydrin Method
9	CLINICAL BIOCHEMISTRY	Dried Blood Spot (DBS)	Pregnancy Associated Plasma Protein A (PAPP-A)	Time Resolved Fluoroimmunoassay
10	CLINICAL BIOCHEMISTRY	Maternal Serum	Beta - Human Chorionic Gonadotropin (Free β -hCG)	Time Resolved Fluoroimmunoassay
11	CLINICAL BIOCHEMISTRY	Maternal Serum	Dimeric Inhibin-A	Time Resolved Fluoroimmunoassay
12	CLINICAL BIOCHEMISTRY	Maternal Serum	Free Beta - Human Chorionic Gonadotropin (Free β -hCG)	Time Resolved Fluoroimmunoassay
13	CLINICAL BIOCHEMISTRY	Maternal Serum	Human Alpha Fetoprotein (hAFP)	Time Resolved Fluoroimmunoassay
14	CLINICAL BIOCHEMISTRY	Maternal Serum	Placental Growth Factor (PlGF)	Time Resolved Fluoroimmunoassay
15	CLINICAL BIOCHEMISTRY	Maternal Serum	Pregnancy Associated Plasma Protein A (PAPP-A)	Time Resolved Fluoroimmunoassay
16	CLINICAL BIOCHEMISTRY	Maternal Serum	Unconjugated Estriol (uE3)	Time Resolved Fluoroimmunoassay
17	CYTOGENETICS	Peripheral Venous Blood, Cord Blood, Products of Conception, Chorionic Villus, Amniotic Fluid, Fetal Blood	Fluorescence in situ hybridization (FISH) - Analysis of Chromosome 13, 18, 21, X & Y	FISH - Fluorescence Microscopy
18	CYTOGENETICS	Peripheral Venous Blood, Cord Blood, Products of Conception, Chorionic Villus, Amniotic Fluid, Fetal Blood	Fluorescence in situ hybridization (FISH) - Analysis of Chromosome 21	FISH - Fluorescence Microscopy
19	CYTOGENETICS	Peripheral Venous Blood, Cord Blood, Products of Conception, Chorionic Villus, Amniotic Fluid, Fetal Blood	Fluorescence in situ hybridization (FISH) - Analysis of Chromosome 22q Microdeletion	FISH - Fluorescence Microscopy



National Accreditation Board for Testing and Calibration Laboratories

SCOPE OF ACCREDITATION

Laboratory Name :

REVVITY HEALTH SCIENCES PRIVATE LIMITED, BLOCK NO. 702-704, 7TH FLOOR, TICEL BIO PARK, PHASE II, CSIR RD, THARAMANI, CHENNAI, TAMIL NADU, INDIA

Accreditation Standard

ISO 15189:2012

Certificate Number

MC-6167

Page No

2 of 3

Validity

04/12/2023 to 22/11/2025

Last Amended on

-

S.No	Discipline / Group	Materials or Products tested	Component, parameter or characteristic tested / Specific Test Performed / Tests or type of tests performed	Test Method Specification against which tests are performed and / or the techniques / equipment used
20	CYTOGENETICS	Peripheral Venous Blood, Cord Blood, Products of Conception, Chorionic Villus, Amniotic Fluid, Fetal Blood	Fluorescence in situ hybridization (FISH) - Analysis of Chromosomes X & Y	FISH - Fluorescence Microscopy
21	CYTOGENETICS	Peripheral Venous Blood, Cord Blood, Products of Conception, Chorionic Villus, Amniotic Fluid, Fetal Blood	Karyotyping (GTG Banding, Quinacrine Banding, C Banding & Ag NOR Banding)	Chromosomal Analysis/Light Microscopy
22	MOLECULAR TESTING	Amniotic Fluid (AF), Chorionic Villi (CVS), DNA, DNA-AF, DNA-CVS, Venous Blood, Product of Conception, Cord Blood, Fetal Blood, Dried Blood Spot, DNA-Venous Blood, DNA - Fetal Blood, DNA - Cord Blood	Sanger Sequencing of single specific genes HBB for thalassemia, CYP21A2, ABCD1, GBA; Single Site sequencing for previously identified variants on NGS testing	Sanger Sequencing
23	MOLECULAR TESTING	Amniotic Fluid (AF), Chorionic Villi (CVS), DNA, DNA-AF, DNA-CVS, Venous Blood, Product of Conception, Cord Blood, Fetal Blood, Dried Blood Spot, DNA-Venous Blood, DNA - Fetal Blood, DNA - Cord Blood	SMA - MLPA	MLPA
24	MOLECULAR TESTING	Amniotic Fluid (AF), Chorionic Villi (CVS), DNA, DNA-AF, DNA-CVS, Venous Blood, Product of Conception (POC), DNA-Venous Blood, DNA - POC	KaryoLite BoBs	Bacs-on-Beads
25	MOLECULAR TESTING	Amniotic Fluid (AF), Chorionic Villi (CVS), DNA, DNA-AF, DNA-CVS, Venous Blood, Product of Conception, Cord Blood, Fetal Blood, Dried Blood Spot, DNA-Venous Blood, DNA - Fetal Blood, DNA - Cord Blood	Prenatal BoBs	Bacs-on-Beads
26	MOLECULAR TESTING	DNA, DNA- CV, DNA-AF, Whole Blood, Product of Conception, Chorionic Villi (CVS), Amniotic Fluid (AF), Venous Blood, Fetal Blood, Cord Blood, Cultured Cell- AF, Cultured Cell- CV, DNA- Fetal Blood, DNA- Product of Conception	Maternal Cell Contamination	Testing using Short Tandem Repeat Markers
27	MOLECULAR TESTING	DNA, DNA- CV, DNA-AF, Whole Blood, Product of Conception, Chorionic Villi (CVS), Amniotic Fluid (AF), Venous Blood, Fetal Blood, Cord Blood, Cultured Cell- AF, Cultured Cell- CV, DNA- Fetal Blood, DNA- Product of Conception	Quantitative Fluorescent - Polymerase Chain Reaction	Testing using Short Tandem Repeat Markers
28	MOLECULAR TESTING	Venous Blood	Non- Invasive Prenatal Testing for Trisomy 13,18 and 21	Rolling Circle Replication



National Accreditation Board for Testing and Calibration Laboratories

SCOPE OF ACCREDITATION

Laboratory Name :

REVVITY HEALTH SCIENCES PRIVATE LIMITED, BLOCK NO. 702-704, 7TH FLOOR, TICEL BIO PARK, PHASE II, CSIR RD, THARAMANI, CHENNAI, TAMIL NADU, INDIA

Accreditation Standard

ISO 15189:2012

Certificate Number

MC-6167

Page No

3 of 3

Validity

04/12/2023 to 22/11/2025

Last Amended on

-

S.No	Discipline / Group	Materials or Products tested	Component, parameter or characteristic tested / Specific Test Performed / Tests or type of tests performed	Test Method Specification against which tests are performed and / or the techniques / equipment used
29	MOLECULAR TESTING	Venous Blood	Y- Chromosome Microdeletion	PCR Amplification of Sequence-Tagged-Sites

