

Date

Reference

2022-12-22

2022/474

Scope of accreditation

Testing laboratory according to SS-EN ISO 15189:2012

Revvity Omics Sweden AB

Sollentuna

Accreditation number

10538

A015057-001

Clinical genetics

<i>Technical area</i>	<i>Parameter</i>	<i>Method</i>	<i>Technique</i>	<i>Material</i>	<i>Field</i>	<i>Note</i>
Molecular genetic diagnostics	Probability for Trisomy 13,18 and 21 in fetus (aneuploidy) NIPT	Inhouse method; SOL-SOP-TEST-01 v. 4	Array analysis	Blood	No	Vanadis
			Array analysis	Plasma	No	Vanadis
	Probability for Trisomy 13,18 and 21 in fetus (aneuploidy) NIPT with fetal sex determination	Inhouse method; SOL-SOP-TEST-01 v. 4	Array analysis	Blood	No	Vanadis
			Array analysis	Plasma	No	Vanadis

Clinical immunology and transfusion medicine

<i>Technical area</i>	<i>Parameter</i>	<i>Method</i>	<i>Technique</i>	<i>Material</i>	<i>Field</i>	<i>Note</i>
Clinical immunology	Placental Growth Factor (PIGF)	Inhouse method; SOL-SOP-INST-07 v. 2	Fluoro immunometric assay	Blood	No	Delfia Xpress
			Fluoro immunometric assay	Serum	No	Delfia Xpress

Changes in the scope of accreditation are in bold.