

Certificate

SKQS - Slovak Association for Quality and Management Systems, Ltd.
Alexandra Rudnaya 23, 010 01 Žilina

SKQS confirms on the basis of the audit that:



MatTek in Vitro Life Science Laboratories, s.r.o.

Mlynské Nivy 73, 821 05 Bratislava
Slovak Republic

scope of application

- development, production, marketing and sales of reconstituted tissues of human cell origin for testing the effects of substances, mixtures and formulations,
- research and development in the area of biological sciences and *in vitro* toxicology,
- informative, non-clinical testing of substances, mixtures and formulations *in vitro* (measurements, analysis and controls),
- theoretical and practical professional consulting in the field of *in vitro* toxicology,
- internal and external trainings in the field of *in vitro* toxicology,
- distribution and sales of glass bottom dishes and plates for scientific and research purposes.

has implemented and maintained the management system
in conformance with the requirements of the international standard

ISO 9001:2015

Quality management system

Certificate No.:	568-3/19
Scope number:	23, 34, 35
Validity date:	01. 06. 2022
Issue date:	02. 06. 2019



Ing. Juraj Kotrbanec
Vedúci certifikačného orgánu SKQS