

## Mismatch Repair Deficiency (dMMR) & Microsatellite Instability (MSI-H) Testing in Endometrial Cancer Webinar

**24 September 2024**

**Goals:** It is the aim of this webinar to present an update about MSI/dMMR analysis in Endometrial Cancer from pathologist's perspective in Europe.

### Webinar Program

Start	Duration	Tuesday, 24 September 2024	Referent
16:00	15 min	Welcome	Discovery
16:15	20 min	Rational for an Algorithm-based Approach	Prof. J. Rüschoff
16:35	20 min	Comparison of Different MSI Testing Methods In Endometrial Cancer	Prof. R. Büttner
16:55	20 min	Quality Issues and Quality Assurance	Prof. F. Penault-Llorca
17:15	45 min	Interactive Case Evaluation and Discussion (8 Cases)	Prof. R. Büttner Prof. F. Penault-Llorca Prof. J. Rüschoff
18:00		End of Webinar	

*We assure that the contents of the webinar are designed to be product- and service-neutral. We assure that potential conflicts of interest of the organizer, the scientific management, and the speakers are disclosed to the participants in a self-declaration (e.g., first slide of presentations, handouts, notices, information in the program, link or download). The event (September 24th, 2024) is sponsored by GlaxoSmithKline with €45,600.00.*