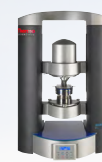


Rheometers & viscometers

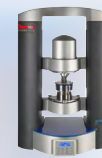
Industry standards compliance

Which rheometer fulfills which industry standard? Entry-level rheometer for absolute viscosity measurements with integrated temperature control. Advanced rheometers for extended material characterization over the largest measuring range.

The following table provides an overview of the compliance of various Thermo Scientific™ HAAKE™ Rheometers with selected industry standards.



| Standard | Industry / sample type | HAAKE Viscotester iQ | HAAKE Viscotester iQ Air | HAAKE MARS iQ | HAAKE MARS iQ Air | HAAKE MARS 40 / 60 | Required accessory / measuring geometry |
|--|---|----------------------|--------------------------|---------------|-------------------|--------------------|---|
| ASTM® D2556 | Adhesives (apparent viscosity) | ● | ● | ● | ● | ● | Viscometers spindles for relative viscosity, universal container holder |
| ASTM D4092 (similar to ISO® 6721-1) | Polymers (Standard terminology for dynamic mechanical properties of plastics) | ✘ | ✘ | ● | ● | ● | Parallel plates or solids clamping tool, parallel plates, upper and lower electrical temperature modules, or temperature chamber |
| ASTM D4440 (equivalent to ISO 6721-10) | Polymers (oscillatory measurements) | ✘ | ✘ | ● | ● | ● | Parallel plates, upper and lower electrical temperature modules or temperature chamber |
| ASTM D4473 | Polymers (oscillatory measurements, curing of thermosetting resins) | ✘ | ✘ | ✘ | ● | ● | Parallel plates, TM-PE-P, PEEK cover, TM-IN-H or TM-EL-H depending on temperature range. Cone and plate only in the case of isothermal tests. Solids clamping tool for composites |



| Standard | Industry / sample type | HAAKE Viscotester iQ | HAAKE Viscotester iQ Air | HAAKE MARS iQ | HAAKE MARS iQ Air | HAAKE MARS 40 / 60 | Required accessory / measuring geometry |
|---------------------------------------|---|----------------------|--------------------------|---------------|-------------------|--------------------|--|
| ASTM D5279 (equivalent to ISO 6721-7) | Polymers (torsional DMA) | ✗ | ✗ | ✗ | ● | ● | Solids clamping tool, temperature chamber |
| ASTM D7271 | Paints & Coatings (oscillatory rheology) | ✗ | ✗ | ✗ | ● | ● | Cone C20 2°, TM-PE-P, PEEK hood |
| DIN® 1342 1-3 | General rheometry | ● | ● | ● | ● | ● | Absolute measuring geometries, temperature control |
| DIN 51810-1 | Lubricating greases | ✗ | ✗ | ✗ | ● | ● | Cone C25 1°, TM-PE-P, PEEK cover |
| DIN 51810-2 | Lubricating greases | ✗ | ✗ | ✗ | ● | ● | Plate P25, TM-PE-P, PEEK cover (tests at +25 °C), TM-IN-H and AC200 thermostat for tests at -40 °C |
| DIN 51810-4 | Lubricating greases | ✗ | ✗ | ✗ | ● | ● | Cone C25 1°, TM-PE-P, PEEK cover |
| DIN 53019-1 | General rheometry (coaxial cylinder measuring geometry) | ● | ● | ● | ● | ● | Coaxial cylinder (with cone at bottom of rotor) |
| DIN 53019-2 | General rheometry (calibration and uncertainty) | ● | ● | ● | ● | ● | Coaxial cylinder (with cone at bottom of rotor) or cone and plate |
| DIN 53019-3 | General rheometry | ● | ● | ● | ● | ● | Absolute measuring geometries |
| DIN 53019-4 | General rheometry (oscillatory rheology) | (●) | ● | ● | ● | ● | Absolute measuring geometries |
| DIN EN 17408 (replaced DIN 54458) | Adhesives | ✗ | ✗ | ✗ | ● | ● | Plate P25, TM-PE-P, PEEK cover |
| DIN EN ISO 3219 1-2 | General rheometry | (●) | ● | ● | ● | ● | Absolute measuring geometries, temperature control |
| DIN EN ISO 6721-1 | Polymers (oscillatory measurements) | ✗ | ✗ | ● | ● | ● | Parallel plates or solids clamping tool, upper and lower electrical temperature modules or temperature chamber |
| ISO 2884 -1 | Paints and varnishes | ● | ● | ● | ● | ● | Cone & Plate measuring geometry |
| ISO 2884-2 | Paints and varnishes | ● | ● | ● | ● | ● | Relative measuring geometries (disc, balls), in combination with 250 ml tin can |

● Oscillatory measurements are an optional feature



| Standard | Industry / sample type | HAAKE Viscotester iQ | HAAKE Viscotester iQ Air | HAAKE MARS iQ | HAAKE MARS iQ Air | HAAKE MARS 40 / 60 | Required accessory / measuring geometry |
|--|--------------------------|----------------------|--------------------------|---------------|-------------------|--------------------|--|
| ISO 6721-7 | Polymers (torsional DMA) | ✗ | ✗ | ✗ | ● | ● | Solids clamping tool, temperature chamber |
| ISO 6721-10 | Polymers (DMA) | ✗ | ✗ | ● | ● | ● | Parallel plates, upper and lower electrical temperature modules or temperature chamber |
| ICA / IOCCC | Chocolate | ● | ● | ● | ● | ● | Coaxial cylindrical CC25 DIN, TM-PE-C, PEEK cover |
| European Pharmacopoeia (Ph. Eur.) 2.2.8 & 2.2.10 | Pharma | ● | ● | ● | ● | ● | Coaxial cylinder geometries |
| United States Pharmacopoeia (USP) 912 | Pharma | ● | ● | ● | ● | ● | Cone & plate and coaxial cylinder geometries |

More information regarding other standards:

DIN 53015, ISO 12058 [product specification FL52253](#): HAAKE Falling Ball Viscometer

Variation of bitumen and asphalt-specific standards (ASSHTO®, ASTM, DIN EN, FGSV AL)

[Data sheet D033](#): Dynamic shear rheometer (DSR) for bitumen and asphalt tests—HAAKE MARS iQ Air Rheometer

 We can help you select the right rheometer.

Please contact us at thermofisher.com/rheometercontactus