

Guide to skilled food rheology

From ingredients to final products

thermo scientific

Next page

Introduction

Feeding the world's rapidly growing population with affordable and sustainable food is a key driver in the food industry.

Developing sustainable food is about ensuring that the world's growing population has not just sufficient food but access to high-quality, nutritious food that is affordable, minimizes environmental impact, and meets customer expectations in terms of taste and texture.

Rheological characterization is used to optimize the mechanical properties or sensory perceptions of food products to meet high customer expectations regarding taste or texture as well as simulate processing conditions. The high flexibility of rheometers, with a broad range of application-specific measuring cells for tribology, texture, and interfacial analysis or even combined optical methods, allows for a comprehensive characterization of products.

With decades of application know-how, our worldwide demonstration labs, and scientists from different disciplines including food rheology, extrusion, and electron microscopy, we can assist you in realizing your specific food application goals. Talk to our experts today and learn what options are available to further your advancement in this innovative arena (contact us).

Table of content

Food rheology requirements

We provide an overview of modern food rheometers for quality control and research & development.

Flow behavior of chocolate melts

Learn about reproducible Quality Control of chocolate according to ICA method 46 with a Thermo Scientific[™] HAAKE[™] Viscotester[™] iQ Rheometer with Peltier and Thermo Scientific[™] HAAKE[™] RheoWin[™] Software.

Simple 2D texture mapping of characteristic food samples from rheological data

Examine a scientific poster presenting 2D diagrams in which different rheological parameters have been plotted against each other to map semi-solid food samples and create landscapes that group them according to their characteristic textural properties.

The potential of tribo-rheometry in food

• Introduction webinar to tribo-rheometry

Dive into the world of tribology and learn how to perform tribo-rheological measurement across different industries with rototational rheometers such as a Thermo Scientific[™] HAAKE[™] MARS[™]iQ Rheometer.

• Tribo-rheometry solution for Thermo Scientific[™] HAAKE[™] Rheometers Discover the tribo-rheometry capabilities of HAAKE rheometers for food applications.

Advantage of hyphenated rheo-spectroscopy and rheo-microscopy for food development using Thermo Scientific[™] HAAKE[™] MARS[™] 40/60 Rheometers

- Cocoa butter crystallization measured with a Rheo-Raman setup
 Simultaneous rheological and Raman spectroscopic measurements enable the examination of the isothermal
 crystallization of cocoa butter.
- Investigating heat-induced gelation of whey protein using simultaneous rheology and FTIR spectroscopy.
 Simultaneous rheological and ETIR spectroscopic measurements allow the gelation process of whey protein

Simultaneous rheological and FTIR spectroscopic measurements allow the gelation process of whey protein to be studied.

Food resource center

Comprehensive food development and testing solutions such as extrusion, spectroscopy, rheology, and more techniques support the development and production of foods. Find more information and support from our scientists.

Food rheology requirements

From production to consumption, rheological properties play an important role during the entire life cycle of liquid or semi-solid food formulations. Rheological measurements with different types of instruments are performed in the food industry on a daily base. Here are some common examples:

- Simple single-point viscosity measurements for a fast batch release in production
- Flow curve or yield point measurements in the quality control laboratory
- Comprehensive rheological investigations such as mouthfeel, swallowing, or chewing behavior for the development of new formulations in the research and development department

Processability, stability, and consistency are the attributes that determine consumer perception and thus the overall acceptance of the final product. They can be investigated with various rheological test protocols. In food production, every stage requires different instrument capabilities.

Modern viscometers and rheometers are available to help navigate daily measurement needs from development to QC.

	Viscotester iQ Rheometer	MARS iQ Rheometer	Viscotester iQ Air Rheometer	MARS iQ Air Rheometer	MARS 40/60 Rheometer		
Rotation mode	 ✓ 	~	 ✓ 	 	 ✓ 	Methods	
Creep and recovery	X	X	 ✓ 	 	 ✓ 		
Oscillation model	~	~	 ✓ 	 	~		
Squeeze, break- penetration tests	X	~	X	 	~		
Tribology	X	~	X	 	~	Accs	
Interfacial rheology	X	Х	X	X	~		
Hyphenated methods	X	X	X	X	~		
Low viscous fluids (beverages)			~	V	~	Samples	
Medium to high viscous fluids	X	X	 ✓ 	V	~		
Semisolids (mayonnaise, mustard)	~	~	~	v	~		
Solids (meat analog, cheese)	X	~	X	V	~		
	Quality control			De fo	Development/ formulation		

Ball bearing models

(+) Back to contents (+) Previous page (+) Next page

Air bearing models

Food rheology requirements

Flow behavior of chocolate melts

Simple 2D texture mapping of characteristic food samples from rheological data The potential of tribo-rheometry in food Hyphenated rheo-spectroscopy and rheo-microscopy in food development

Flow behavior of chocolate melts

Viscosity and yield stress hold outsized importance when making chocolate

The flow behavior of molten chocolate is a crucial property for many reasons. During production, the various transportation, filling, dipping, coating, or dosing steps depend on a precisely defined viscosity and yield stress. The finished chocolate's properties, such as its surface's appearance or mouthfeel, are also directly related to the viscosity behavior of the chocolate.

In this application note, the rheological properties of two different chocolate samples were measured according to ICA Method 46 with a HAAKE Viscotester iQ Rheometer with Peltier and HAAKE RheoWin Software.

Read the application note





(+) Back to contents (+) Previous page (+) Next page

Food rheology requirements

Flow behavior of chocolate melts

Simple 2D texture mapping of characteristic food samples from rheological data The potential of tribo-rheometry in food

Hyphenated rheo-spectroscopy and rheo-microscopy in food development

Simple 2D texture mapping of characteristic food samples from rheological data

Categorize thickened food products using distinctive textural attributesin new end products

Overview

Rheological testing is a powerful and well-established tool for mechanically characterizing a range of food samples, including liquids, semisolids, and solids. However, relying solely on a single number, such as the apparent viscosity at a specific shear rate, proves inadequate for fully describing a material's textural properties. A more comprehensive investigation is necessary here, especially if different foods are to be grouped and classified according to their rheological behavior.

This scientific poster examines the results of three rheological tests conducted on nine different semi-solid food products. Taking the right parameters and plotting them against each other in a 2D diagram can help to identify similarities and better understand structure-property relationships of foods.

View the poster





(Back to contents (Previous page (Next page

Food rheology requirements

Flow behavior of chocolate melts Simple 2D texture mapping of characteristic food samples from rheological data

The potential of tribo-rheometry in food Hyphenated rheo-spectroscopy and rheo-microscopy in food development

The potential of tribo-rheometry in food

Know that your new innovative food will have customer acceptance

Tribo-rheometry is a valuable technique used in the food industry to assess the rheological properties of food products. Measuring the frictional forces generated during the mastication and swallowing of food materials helps in understanding their texture, stability, and sensory attributes. This information is crucial for optimizing food formulations, enhancing product quality, and improving consumer satisfaction.

Webinar

Introduction to tribo-rheometry (18 minutes)

Dive into the world of tribology and learn how to perform tribo-rheometrical measurements on rotational rheometers across industries.

Video

Tribo-rheometry solution for Thermo Scientific HAAKE Rheometers

Discover the tribo-rheometry capabilities of HAAKE Rheometers for food applications.

(Watch the webinar





Watch the video





Food rheology requirements

Flow behavior of chocolate melts Simple 2D texture mapping of characteristic food samples from rheological data The potential of tribo-rheometry in food

Hyphenated rheo-spectroscopy and rheo-microscopy in food development

Advantages of hyphenated rheo-spectroscopy and rheo-microscopy in food development See more with multimodal analytical techniques

The combination of two complementary measurement techniques at the same time can reveal more information about a food's properties than either one of them individually. Explore these two application notes to see examples.

Application note

Cocoa butter crystallization measured with a Rheo-Raman setup Simultaneous rheology and Raman spectroscopy measurements enable the examination of the isothermal crystallization of cocoa butter.

Application note

Investigating heat-induced gelation of whey protein using simultaneous rheology and FTIR spectroscopy Simultaneous rheological and FTIR spectroscopic measurements allow the gelation process of whey protein to be studied.

Read the application note





Read the application note

Investigating cocoa butter crystallization unvestigating cocoa putter crystallication using meology and Raman spectroscopy

Investigating heat-induced gelation of whey protein using simultaneous rheology and FTIR spectroscopy

thermo scientific

(Back to contents (Previous page (→) Next page

Food rheology requirements

Flow behavior of chocolate melts Simple 2D texture mapping of characteristic food samples from rheological data

The potential of tribo-rheometry in food Hyphenated rheo-spectroscopy and rheo-microscopy in food development

Thermo Fisher

Food resource center

Thermo Fisher



What rheology can do for food From rheology to mouthfeel

Do you know that rheology plays a crucial role in determining the sensory attributes, stability, and processing characteristics of various food products.

By studying the flow behavior and texture of food, researchers can design formulations and processing conditions to achieve desired sensory attributes and ensure product stability.

Watch the webinar

thermo scientific



thermo scientific

Application laboratories

Our fully equipped application laboratories are in constant demand for testing customer samples and developing and optimizing pioneering applications. We provide a broad range of product and application solutions, and our team of application scientist and interdisciplinary technique specialists is on hand to answer your questions (contact us).

Register for application and product information at <u>thermofisher.com/specoptin</u> to gain access to the latest resources to accelerate your research and improve laboratory productivity.

Seminars and trainings

Comprehensive training programs, in-house seminars, and practical rheology and extrusion courses in various locations around the world. We support our customers with comprehensive on-demand training (webinars, videos, and application notes) from our experts, and benefit from our interdisciplinary knowledge in food science and technologies.

Interested in food sustainability?

<u>Read our compendium</u>, food extrusion and rheology in alternative protein formulations and valorization of food side-streams.

Food rheology: thermofisher.com/foodrheology Food extrusion: thermofisher.com/plantbasedmeat

For research use only. Not for use in diagnostic procedures. For current certifications, visit thermofisher.com/certifications.
© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. PPA-C01046 07/24 M