

# Thermal Analysis Method Development

Authors: Andrew W. Salamon, Dr. Peng Ye PhD., Dr. Kevin Menard PhD.

# INTRODUCTION

This poster guides the new thermal analyst through the most demanding areas of analytical science. The difficulty of defining the appropriate thermal analytical technique and how to develop a thermal method for that technique are simplified.

### Method Development Steps to follow

- . Step 1 -- Ask a series of questions.
- · Step 2 -- Identify the Thermal Analysis Techniques to use
- \* Step 3 -- If applicable, select the appropriate ASTM method to use
- Step 4 -- Enlist expert knowledge to assist and interpret data.

## Step 1 -- Ask a Series of Questions.

This is the ground work for the best thermal technique and method.

#### Ouestion 1 -- What is the physical state of the material?

- Is the material state a powder?
- Is the material state a pellet?
- Is the material state a paste?
- Is the material state a foam?
- Is the material state a thin film?
- Is the material state a fiber? Is the material state a liquid?
- Is the material state a rectangular bar?
- Is the material state an irregular shape?
- Is the material state a tube?
- Is the material state a gel? Is the material state a cylinder?

These questions begin to identify what thermal techniques and accessories should be used.

### Question 2 -- What are the material's physical properties to be analyzed by Thermal Analysis?

Do they include

Soften points?

Effects of plasticizer?

Glass transition temperate?

Melting point? Heat of fusion

Specific Heat?

Purity?

Oxidative stability

Rates of Reaction (Kinetics) Curing?

Cross linking?

Degree of cure?

Decomposition? Optimum Crystallization temperature?

Pyrolysis?

Weight percent filler?

Weight loss vs. time? Weight loss vs. temperature?

Loss of water?

Loss of solvent?

Loss of plasticizer? Impact properties?

Viscoelastic behavior?

Mechanical behavior?

Long term behavior?

Modulus vs. temperature? Damping vs. temperature?

Chain branching?

Molecular weight?

Molecular weight distribution?

Liquidus or Solidus temperatures?

# Question 3 -- What type of information is required?

Quantitative' Qualitative?

A mixture of both?

#### Ouestion 4 -- What analytical accuracy and precision is required?

Accuracy is the extent to which the results of a measurement approach the

Precision is measure of the range of values of a set of measurements

Example: When five darts are thrown at a bull's eye hit the bull's eye, then the accuracy is said to be good. The fact that all five darts are close to each other indicates good precision.

Question 5 -- Are the events of interest minor or major?

Question 6 -- What do you already know about the material or sample?

Question 7 -- What are the important properties of the material?

Question 8 -- Is the material a single component or a complex mixture?

Question 9 -- What is the material's history?

Question 10 -- What is the material's future?

Question 11 -- Is there a limitation on sample size?

Question 12 -- How many samples must be run?

Ouestion 13 -- What is the required analysis turn-around time?

Ouestion 14 -- Are there any safety hazards to be concerned about?

# Step 2 – Use the applications listed below to help identify appropriate Thermal Analysis techniques for your investigation. Many times a combination of techniques are required to solve the problem.

Differential Scanning Calorimetry (DSC) is the workhorse of thermal analysis laboratory. DSC can examine materials between -180C to 750C.

**Definition:** DSC is a technique upon which the heat flow to or from a sample specimen is measured as a function of temperature or time as the sample specimen is subjected to a controlled temperature program in a controlled atmosphere.

## Some DSC applications are listed below:

- \*Identify the softening point of a material (glass transition)
- •Compare additive effects on a material
- ·Identify the glass transition temperature.
- \*Identify a material's minimum process temperature.
- \*Identify the amount of energy to melt a material.
- ·Quantify a material's specific heat.
- \*Perform oxidative stability testing (OIT).
- \*Characterize a material as it cures under UV light.
- . Optimize the crystallization temperature upon cooling.
- \*Study polymorphic stability in pharmaceuticals

#### Some more DSC applications are listed below:

- Study amorphous content.
- . Understand the reaction kinetics of a thermoset material as it cures.
- . Compare the degree of cure of one material to another
- \*Examine thermal events that are very week or masked by other material characteristics by employing very fast scanning rates (Hyper-DSC ®).

# Thermogravimetric Analysis (TGA) examines materials between ambient and

Definition: TGA is a technique in which the mass of a substance is monitored as a function of temperature or time as it is subjected to a controlled temperature program in a controlled atmosphere.

#### Some TGA applications are listed below:

- ·Identify the filler content of a material by weight percent.
- ·Identify the ash content of a material by weight percent.
- \*Characterize a material's weight loss within a certain temperature range.
- \*Characterize a material's weight loss vs. time at a given temperature.
- Ouantify a material's loss of water, solvent, or plasticizer within a certain temperature range.
- ·Examine flame retardant properties of a material.
- ·Examine the combustion properties of a material.
- ·Perform oxidative stability testing

# Dynamic Mechanical Analysis (DMA) examines materials between

**Definition:** DMA is a technique in which a substance while under an oscillating load is measured as a function of temperature or time as the substance is subjected to a controlled temperature program in a controlled atmosphere.

#### Some DMA applications are listed below:

- ·Quantify impact properties of a material.
- . Examine the Viscoelastic behavior of a material.
- ·Examine a material's mechanical behavior.
- •Examine a material's long term behavior while under load.
- ·Ability to extrapolate material behavior on a time basis.
- ·Identify a material's modulus vs. temperature.
- ·Examine a material's damping qualities vs. temperature.
- ·Examine the effects of temperature on molecular chain branching.
- ·Compare a material's molecular weight and weight distribution. ·Examine additive effects on a material's mechanical properties.
- \*Material -humidity studies from 10% RH to 80% RH

# Step 3 - Research ASTM methods. There may be a traditional ASTM method that is suited for your investigation.

•There are more than 25 ASTM methods for each thermal technique

•There are several ASTM methods that can utilize more than one technique of thermal

# Step 4 - If required, seek out expert knowledge to review your data, suggest additional testing, and recommend possible solutions to your problem.

#### For expert knowledge look to:

- •Local Universities
- •Consultants
- •Instrument manufacturers