

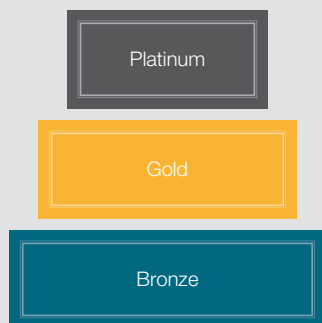
Customer Tiered Support Model

A model to ensure our customers success

What is a customer tiered support model?

Thermo Fisher™ Platform for Science™ software utilizes a tiered support model to ensure that we are offering support that is tailored to meet the needs of our customers in production. As our customers' needs grow, they will move up to the next tier.

What are the tiers?



Some examples of how a tier is defined

- Multi-site or Single-site expansion
- Number of registered users or annual financial commitment
- Business factors such as new market configurations

Key benefits by tier

Platinum	Gold	Bronze
<ul style="list-style-type: none">• Senior Customer Success Manager to monitor user adoption, train on configuration and product capabilities, triage configuration related issues, evaluate expansion opportunities, and maintain communication at the user/manager level• Senior Account Manager to manage renewals and subscription changes• Executive Business Review (EBR) to verify system adoption and define future plans for expansion• Thermo Fisher Stakeholder to establish and maintain routine communication with customer stakeholder• Customer Support Team to assist on all questions and facilitate system upgrades	<ul style="list-style-type: none">• Customer Success Manager to monitor user adoption, train on configuration and product capability, triage configuration related issues, evaluate expansion opportunities, and maintain communication at the user/manager level• Account Manager to manage renewals and subscription changes• Thermo Fisher Stakeholder to establish and maintain routine communication with customer stakeholder, and provide strategic direction for future phases of expansion• Customer Support Team to assist on all questions and facilitate system upgrades	<ul style="list-style-type: none">• Tier 3 Specialist to maintain communication at the user/manager level• Customer Support Team to assist on all questions and facilitate system upgrades• Thermo Fisher Stakeholder available as needed

Find out more at coreinformatics.com