

Supplier Diversity Program Frequently Asked Questions

1. Which diverse classifications does Thermo Fisher Scientific accept in the Supplier Diversity Program?

Our Supplier Diversity Program provides qualified certified small and diverse businesses with a pathway to become an official provider of goods or services for Thermo Fisher Scientific.

Suppliers that qualify for this program include:

- Minority Owned Business
- Women Owned Business
- LGBT Owned Business
- Veteran Owned Business
- Service-Disabled Veteran Owned Business
- Disabled Owned Business
- Small Businesses registered through the Small Business Administration (SBA)

2. Which certifying agencies does Thermo Fisher Scientific accepts?

Supplier Diversity Program accepts all State, and local government Supplier Diversity certifications. In addition, we accept most major national/regional councils such as:

- National Minority Supplier Development Council (NMSDC) www.nmsdc.org
- Women's Business Enterprise National Council (WBENC) www.wbenc.org
- National Veteran-owned Business Association (NaVOBA) www.navoba.com
- National Veteran Business Development Council (NVBDC) www.nvbdc.org
- National Gay and Lesbian Chamber of Commerce (NGLCC) www.nglcc.org
- Disability:IN www.DisabilityIn.org
- Diversity Alliance for Science www.diversityallianceforscience.com

3. How does a supplier do business with Thermo Fisher Scientific through its Supplier Diversity Program?

To be considered for business through the Supplier Diversity Program, your company must meet the certification and registration requirements as one of the supplier classifications shown in the previous answer. If your company has met our certification requirements, we encourage you to register with our program through [VIVA STARS](#)



4. What happens after I register with the Supplier Diversity Program?

After you've completed your registration, your information will reside in our database that is provided to sourcing/procurement staff within Thermo Fisher Scientific. If there's a sourcing opportunity, our buyers and sourcing staff will access this database to search for suppliers based on goods and services, and geographical location information.

5. Who do I contact if I have questions about my STARS (VIVA) registration or technical questions?

- For any technical issues please reach out via email to support@starssmp.com or via phone at (877) 557-5258.

6. What are other actions or next steps I can take to be more proactive?

Thermo Fisher Scientific company buyers annually attend outreach events and supplier trade shows across the United States. These events are listed on the [Outreach page of the website](#).

7. If I have questions about the Supplier Diversity Program?

If after reading the Supplier Diversity Program section of the website you have additional questions please send an email to supplierdiversity@thermofisher.com