

PerkinElmer Financial Services
Powered by Wells Fargo

Complete Financing Solutions



PerkinElmer Financial Services understands the environment you are competing in and the power a new equipment solution can play in increasing your productivity and competitiveness. That's why we provide fast, simple and reliable financing solutions that help you achieve your business goals.

Unleash the power of financing

Whatever objectives are driving your business, financing can help you capitalize on opportunities, while staying on budget.

Financing can offer:

- Predictable and affordable payments
- Reduced up-front costs and preserved working capital
- Customized structures and tailored payment solutions
- Easy equipment and technology upgrades
- Potential tax and accounting advantages

Our Financing Products

Leases, such as the Fair Market Value Lease, can offer end of term flexibility with a clear path for your next equipment upgrade.

Loans can provide low monthly payments while maintaining the benefits of ownership and associated tax benefits. These options are ideal for customers with long-life equipment needs or businesses who want to structure monthly expenses to match their cash flow.

PerkinElmer Financial Services also offers specialty products designed to meet the unique needs of your industry. Payments can be structured in many ways including: monthly, quarterly, deferred, step, seasonal and extended terms.

Tools, resources & expertise to help you build a stronger business

Speed

PerkinElmer Financial Services knows you are busy and financing shouldn't slow you down. We offer a fast, efficient credit process — often with a credit decision within minutes — to get you funded and out the door as quickly as possible.

Convenience

Financing from PerkinElmer Financial Services is tailored to meet your needs. Here's what that could look like for you:

Financing can offer:

- Consolidate your entire purchase — from equipment to software to installation - into a single convenient payment.
- Avoid the large upfront costs and risks associated with ownership with our options to match a lease term with the length of a job contract or new product development.
- Structure your payments according to cash flow, with options like deferred or stepped payments.
- Structure your lease to take advantage of the best years of the equipment with the flexibility to upgrade at the end of the lease terms.
- Take the hassle out of paying your bills with our easy-to-use online tools.

Flexibility

We offer products and services customized to your cash flow needs. Regardless of whether you are looking to optimize depreciation, lower monthly payments or structure them to meet your business needs, our products and solutions can help.

Expertise

Our people have been doing this for decades; we know how to match our financing solutions to your business needs. When you work with PerkinElmer Financial Services, you're working with financing experts who are committed to adding real value and helping your business grow.

*Call PerkinElmer today to
learn what we can do for
your business.*