

WAVEcare

Smooth operations

WAVEcare PROGRAMS

Carefree or basic support for high-quality kinetic data and unparalleled performance

THE FIRST 12 MONTHS

Care for the first year of your WAVEsystem is on us, with a warranty period that covers travel, labor and parts costs for one (1) on-site preventive maintenance (PM) visit to be carried out within 12 months. Travel, labor and parts costs for unscheduled repairs are also included.

Maintenance is carried out by our qualified service professionals.

WHY PREVENTIVE MAINTENANCE (PM) VISITS?

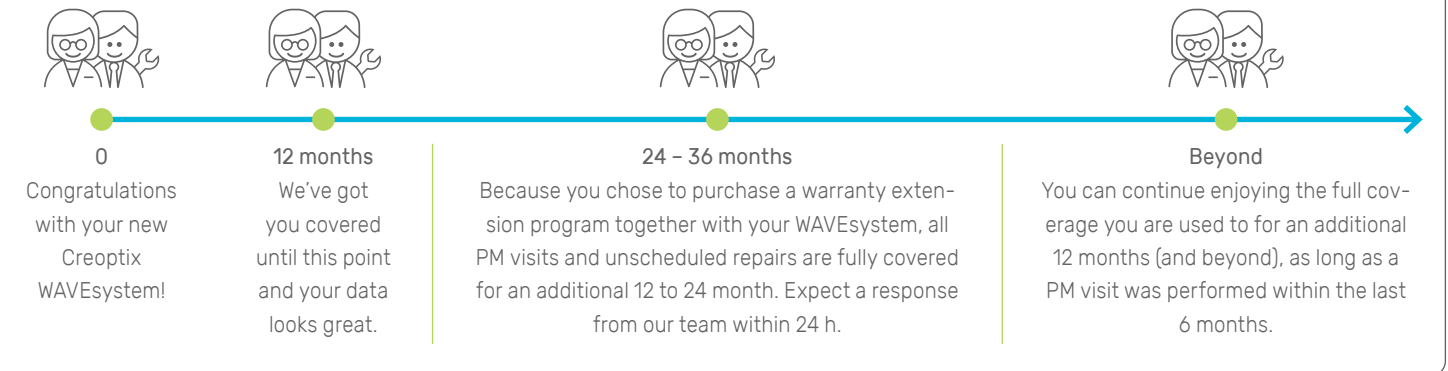
Many fluidic components undergo wear and tear over time. PM visits are crucial for the continuous production of high-quality data, and our team of engineers strongly recommend scheduling one every six months. During a PM visit, a certified member of our team of engineers will run a full check of the WAVEsystem to assess its performance, including:

- Replacement of various fluidic components
- Calibration of the device
- Update of both firmware and software (applicable only for carefree programs)
- A system function test (SFT), including an official report

The different WAVEcare programs ensure the continuous, high performance operation of your WAVEsystem. Choose the program that fits your needs.

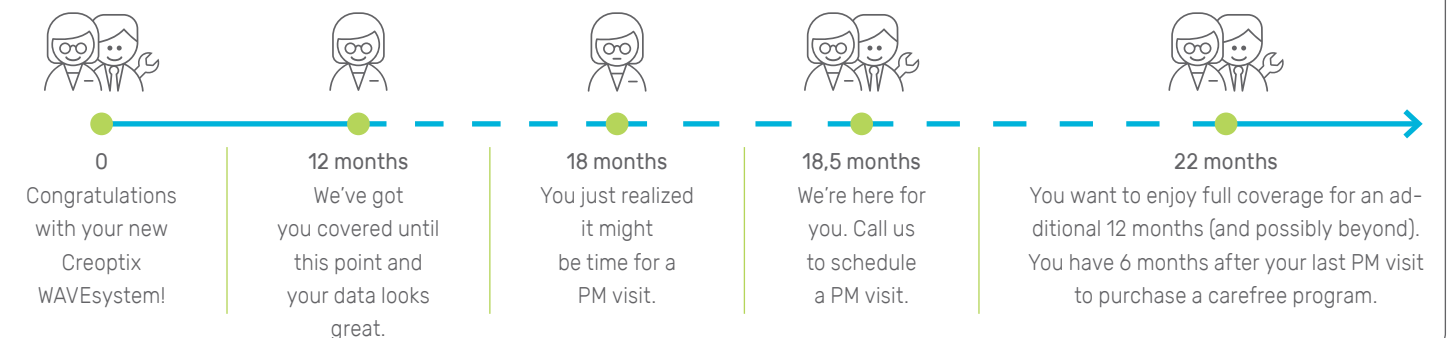
CAREFREE PROGRAM: for the regular user

Know what you are up to and plan your maintenance budget ahead of time



BASIC PROGRAM: for the occasional user

Call us when you need us the most for a PM visit or a repair



ORDERING INFORMATION

Choose any of the WAVEcare programs below to complement the initial 12 month warranty period included with your new WAVEsystem

CARE-FREE: ONGOING SUPPORT FOR THE REGULAR USER

Response time: up to 24 hours

ORDER NO.	PM visits		Travel, labor and parts costs for unscheduled repairs	Software upgrades	Time of purchase	Validity
	Number included	Travel, labor and parts costs				
[WARR-3Y] Warranty extension to 36 months	5	✓	✓	✓	Before installation	36 months
[WARR-2Y] Warranty extension to 24 months	3	✓	✓	✓	Before installation	24 months
[FULL-1Y] Full coverage for 12 months	2	✓	✓	✓	Within 6 months after a PM visit	12 months

BASIC: A LA CARTE SUPPORT FOR THE OCCASIONAL USER

Response time: up to 5 working days

ORDER NO.	PM visits		Travel, labor and parts costs for unscheduled repairs	Software upgrades	Time of purchase	Validity
	Number included	Travel, labor and parts costs				
[MAINT-1PM] Maintenance contract - 1 PM	1	✓	✗	✗	Anytime	12 months
[MAINT-2PM] Maintenance contract - 2 PMs	2	✓	✗	✗	Anytime	12 months
[VISIT-PM] Preventive maintenance visit	1	✓	✗	✗	Anytime	4 weeks

EXTRA'S

ORDER NO.	Description
[LABOR-1h] Labor hourly rate	On-site labor from one of our service team members.
[RELOCATE] Relocation service	Relocation service for the WAVEsystem, including system function test (SFT).
[REMSUP-1h] Remote support hourly rate	Remote support from one of our service team members.
[TRAVEL-1h] Travel hourly rate	Travel costs for one of our service team members.

WAVEcare programs apply to all WAVEsystems. For special requests, contact us through support@creoptix.com

Place your order through orders@creoptix.com

WE FOCUS ON SMOOTH OPERATIONS, YOU FOCUS ON DELIVERING RESULTS

To stay up-to-date with our latest developments,
visit www.creoptix.com

To make the most out of your WAVEsystem,
contact our team at support@creoptix.com



LABEL-FREE DATA
LIKE YOU'VE NEVER SEEN BEFORE



Creoptix AG | Zugerstrasse 76 | 8820 Wädenswil | Switzerland (HQ)
Creoptix Inc | c/o The Brighton Company | 111 South Bedford Street, Ste. 108 | MA-01803 Burlington | USA
www.creoptix.com | +41 44 533 26 60

4013BR 01/2020 2

©2020 Creoptix Sensors / Creoptix, WAVEchips, WAVEcore and waveRAPID are trademarks of CreoptixAG. All goods and support services sold by Creoptix are subject to our terms and conditions. Please visit our website or contact your local Creoptix representative for the most up-to-date information.

KEEPING KINETICS REAL