

How to read your Dosimetry Report

Periodic reports are now just that. If you are looking for information on a "June dosimeter", you will find it on your June report, regardless of the dosimeter return date.

When returned with your dosimeters, the control dosimeter lists values in the calculation of your assigned dose. YTD and Lifetime are not tracked as these values in themselves are not actual dose.

ND - indicates that your dosimeter was processed, but the dose is below our minimum reportable dose of 10 millirem.
NR - Indicates that the dosimeter does not monitor dose in that category.

Highlighted areas contain your assigned/effective dose for this report. It is representative of either a single dosimeter dose, the calculated dose of multiple dosimeters listed immediately underneath, or an Effective Dose Equivalent (EDE) calculation.

Users working in multiple groups should be tracked across all groups to ensure compliance to yearly exposure limits.

ALARA Notifications can now be found at the end of each report. If you have a specific ALARA limit you would like to have set for you, please contact a Customer Care representative. We will be happy to make that adjustment for you.

Each report generated has a version number. Always viewing the version with the highest number ensures you are looking at the most timely, accurate report.

ACCOUNT 12345 GROUP 2 WEAR PERIODS ENDING JUNE 2018 VER 2 PUBLISHED 07/10/2018 REPORT NO. 20012514
QA CASE ACCOUNT / QA GROUP 2

MONTHLY DOSIMETRY REPORT

PIN	NAME	CALCULATION	DOSIMETER TYPE	LOC	WEAR PERIOD	NOTES	CURRENT DOSE (MILLREM)			DOSE THIS YEAR (MILLREM)			LIFETIME DOSE (MILLREM)			
							NEU	DEEP	EYE	SHALLOW	DEEP	EYE	SHALLOW	DEEP	EYE	SHALLOW
0001014	CONTROL	Control	82	XC	06/01/2016-06/30/2018	Control	NR	ND	ND	ND	ND	ND	ND	ND	ND	ND
0001009	ANDERSON ALAN	Standard	82	CH	06/01/2016-06/30/2018	URN	Body dose				ND	ND	ND	ND	ND	ND
0001011	DAWSON HARRY	Standard	82	CH	06/01/2016-06/30/2018		Body dose	ND	ND	ND	ND	ND	ND	10	ND	ND
		Standard	82	CH	06/01/2016-06/30/2018		Extremity dose				ND		ND			ND
0001021	GONZALES SUSAN	Standard	05	RU	06/01/2016-06/30/2018		Body dose	125	125	125	125	125	125	125	125	125
0001007	JOHNSON BOB	Standard	82	CH	06/01/2016-06/30/2018		Body dose	NR	ND	ND	ND	11	11	11	17	17
0001008	LEE LORETTA	Standard	82	CH	06/01/2016-06/30/2018		Body dose	NR	ND	ND	ND	ND	ND	ND	SEE MULTIGROUP	
0001002	RICHARDS CECIL	Standard	82	CH	06/01/2016-06/30/2018		Body dose	NR	ND	ND	ND	ND	ND	ND	SEE MULTIGROUP	
		Webster 1	82	CL	06/01/2016-06/30/2018		Body dose	NR	10	10	10	ND	10	10	ND	10
0001001	WILLIAMS WILMA	Webster 2	82	CL	06/01/2016-06/30/2018		Body dose	NR	10	10	10	24	30	30	24	30
		Webster 2	82	WS	06/01/2016-06/30/2018		Body dose	NR	10	10	10	ND	ND	ND	ND	ND

Explanation of Notes
UNR UnReturned Badge

MultiGroup
The following wearers have been monitored in multiple wear groups. This section summarizes the results from all wear groups for each wearer.

PIN	NAME	CALCULATION	DOSIMETER TYPE	LOC	WEAR PERIOD	NOTES	DOSE THIS YEAR (MILLREM)			LIFETIME DOSE (MILLREM)		
							NEU	DEEP	EYE	SHALLOW	DEEP	EYE
0001008	LEE LORETTA	Body dose					10	10	10	10	10	10

Notes will be given an alpha or numeric code which is referenced at the end of the report for explanation. An example would be this "UNR" used as unreturned

ALARA Notifications
The following wearers have received a dose that exceeds an ALARA threshold. See the threshold values on the final page of this report.

PIN	NAME	THRESHOLD	CURRENT DOSE (MILLREM)			
			DEEP	EYE	SHALLOW	
0001021	GONZALES SUSAN	Level I	Body dose	125	125	125

Explanation of Notes
UNR UnReturned Badge

ACCOUNT	GROUP	WEAR PERIODS ENDING	VER	PUBLISHED	REPORT NO.											
12345	2	JUNE 2018	2	07/10/2018	20012514											
QA CASE ACCOUNT / QA GROUP 2																
PIN	NAME	CALCULATION	DOSIMETER TYPE	LOC	WEAR PERIOD	NOTES	NEU	DEEP	EYE	SHALLOW	DEEP	EYE	SHALLOW	DEEP	EYE	SHALLOW
1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17

- Assigned wearer identification number.
- Employee name, area monitor location, or other reference designation.
- Calculation methodology used for assigned dose computation.
- Dosimeter Type:
82 XBGN
83 XBGN + Track Etch
82W Wrist TLD XBGN
82E Environmental TLD XBGN
83E Environmental TLD XBGN + Track Etch
05 Extremity ring
- Location code - where on the body the dosimeter is worn:
XC Control FM Fetal Monitor
CH Chest EV Environmental or Area Monitor
CL Collar RU Right Hand or Upper Extremity
WS Waist LU Left Hand or Upper Extremity
- The wear period the dosimeter monitored.
- Notes pertaining to the specific dosimeter.
- Neutron component of Deep Dose.
- Deep Dose Equivalent applies to external whole body exposure and is the dose equivalent at a tissue depth of 1 centimeter (1000 mg/cm2). Dose in millirem is reported for photon energies from approximately 10 keV to MeV. Neutron dose is included if applicable.
- Eye Dose Equivalent applies to the external exposure to the lens of the eye and is the dose equivalent at a tissue depth of 0.3 centimeters (300 mg/cm2). It includes dose in millirem for beta particles and photons. Neutron dose is included if applicable.
- Shallow Dose Equivalent applies to the external exposure of the skin or an extremity and is the dose equivalent at the tissue depth of 0.007 centimeters (7 mg/cm2) averaged over an area of 1 square centimeter. It includes dose in millirem for beta particles and photons. Extremity doses are reported in this column based on 662 keV photons unless other energy or radiation source information is available. Neutron dose is included if applicable.
- Year to Date Cumulative Deep Dose. Neutron dose is included if applicable.
- Year to Date Cumulative Eye Dose. Neutron dose is included if applicable.
- Year to Date Cumulative Shallow Dose. Neutron dose is included if applicable.
- Lifetime Cumulative Deep Dose. Neutron dose is included if applicable.
- Lifetime Cumulative Eye Dose. Neutron dose is included if applicable.
- Lifetime Cumulative Shallow Dose. Neutron dose is included if applicable.

Explanation of Dosimetry Report